

GenSmart™ 2.0 Online Ordering

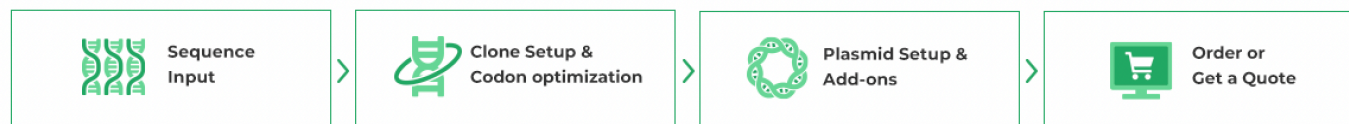
Quick User Guide

Order Gene Synthesis

The GenSmart™ 2.0 online ordering platform is an integrated online order center for all your signature molecular biology services. With this top-notch platform, ordering has never been easier.

Experience a seamless journey as you input, edit, and optimize your gene sequences, while conveniently ordering gene synthesis, cloning and plasmid DNA preparation services tailored to your specific needs. The platform features the best-in-class GenSmart™ codon optimization tool, allowing you to optimize sequences with a single click. Discover the best services at the most competitive prices with GenSmart™ 2.0 now.

Workflow Overview



Order Gene Synthesis via GenSmart™ 2.0 with Ease

- ✓ Utilize the best-in-class codon optimization algorithm to optimize your sequence
- ✓ Select the most suitable gene synthesis services for your applications
- ✓ Receive estimated pricing and turnaround time
- ✓ Get an instant quote and place your order

1. Sequence Input

- Select and click on “Gene Synthesis” tab. Then input your project name and select your downstream application from the drop-down menu. Click “Start New Project” to start gene synthesis online ordering (see **Figure 1**).
- Input your nucleotide and/or amino acid sequences by copying and pasting the sequences into the sequence table manually (see **Figure 2**).
- Or upload sequence files to input your sequences. GenSmart™ 2.0 accepts Excel, CSV/TSV, FASTA and other commonly-used types of files for uploading sequences, and you can find the file templates on the bottom of the page (see **Figure 2**).
- Once sequence input is done, click “Continue” to proceed to the basic setting page for cloning settings and codon optimization.

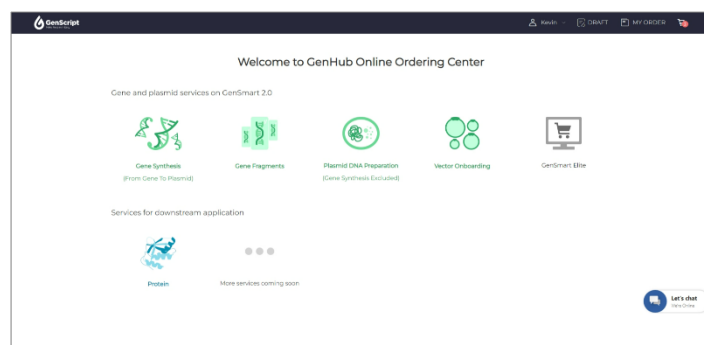


Figure 1: GenSmart™ 2.0 online ordering platform. Order and quote gene synthesis and plasmid DNA preparation services in only 4 simple steps, as well as connect and get support from our Ph.D. level experts.

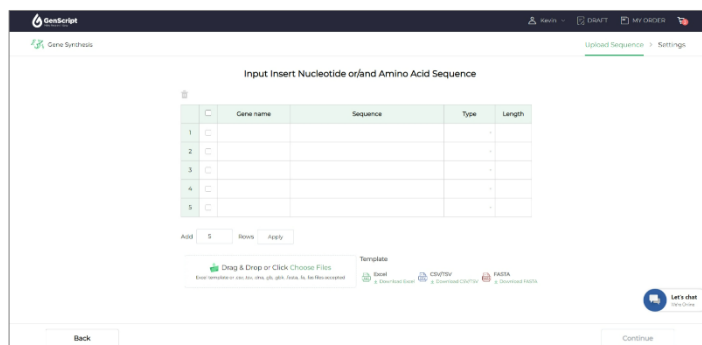


Figure 2: Sequence input. Input nucleotide and/or amino acid sequences by uploading sequence files or manually input. Drag & Drop or choose files to upload the sequence files. The templates of Excel, CSV/TSV, and FASTA files are available on the bottom of the page.

2. Clone Setup and Condon Optimization

- GenSmart™ 2.0 automatically analyzes the synthesis complexity, checks for errors in your sequences, and fills in default options for vectors, 5' and 3' restriction enzymes. Our best-matching program will also select the optimal speed for your sequences and calculate the corresponding price and production time. If everything looks good, click "Continue" to proceed.
- You can also customize the cloning settings. Choose a clone vector, or an expression vector, or a previously archived vector for your sequences. You can apply the chosen vector to one sequence or select "Batch Change Vector" to apply them to all selected sequences (see **Figure 3**).
- Once the vector is applied, choose the cloning method for your sequences from restriction enzyme cloning or seamless cloning in the "5' RE" and "3' RE" columns on the sequence table.

⚠ If you opt for the seamless cloning method, remember to add homology arms to the 5' and 3' ends of your sequences. You can select the corresponding sequences, and then click "Add Flanking" button to add homologous arms to all select sequences (see **Figure 4**). Alternatively, you can click on the specific sequence to add homologous arms in the sequence editing page. For restriction enzyme cloning, choose the restriction enzymes in the 5' and 3' RE columns, and the corresponding sequences will be automatically added to both ends of your sequences.
- Check the "Status" column for the synthesis complexity and errors of your sequences (see **Figure 5**). You can check the complex regions of the complex sequence via the complex icon. For more details, you can click on the complex sequence and check the complex regions in the sequence editing page (see **Figure 6**). We highly recommend performing codon optimization using our best-in-class algorithm to lower synthesis complexity and improve downstream application performance.

⚠ Look out for red error signs and yellow warning signs, as they indicate issues that need to be addressed before processing your order. Click on the signs to view the details.
- Click on the sequence to edit your sequence in the sequence editing page (see **Figure 6**).
- Click "Codon Optimization" to optimize your sequence using best-in-class codon optimization algorithm (see **Figure 6**). Select the region of your sequence that requires codon optimization, the host organism, and the restriction sites to keep and/or avoid and then click "Continue" to start the codon optimization (see **Figure 7**). After successful codon optimizations, you can continue with your original sequence or optimized sequence. You can also download the codon optimization report by clicking on the optimization report icon (see **Figure 8**). Alternatively, you can select on the sequences requiring codon optimization and then click "Batch Optimization" on the basic setting page to preform batch codon optimization (see **Figure 9**).
- GenSmart™ 2.0 will choose the optimal service speed for your sequences. However, you can manually select your preferred speed from FLASH, Premium, Basic, or Rocket. The price and production time will adjust accordingly afterwards (see **Figure 3**).

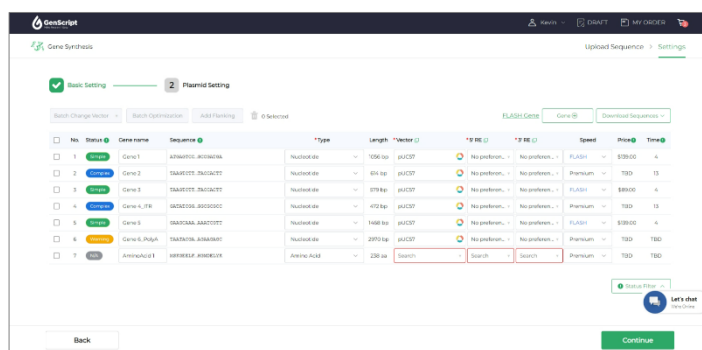


Figure 3: Basic setting page. After the input of the sequences, GenSmart™ 2.0 automatically analyzes the sequences and fills the cloning parameters. You can directly click “Continue” to finish the cloning settings.

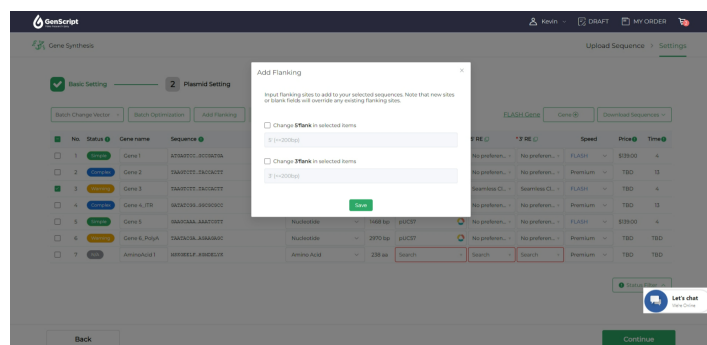


Figure 4: Add homology arms for seamless cloning. Select the corresponding sequences, and then click “Add Flanking” button to add homologous arms.

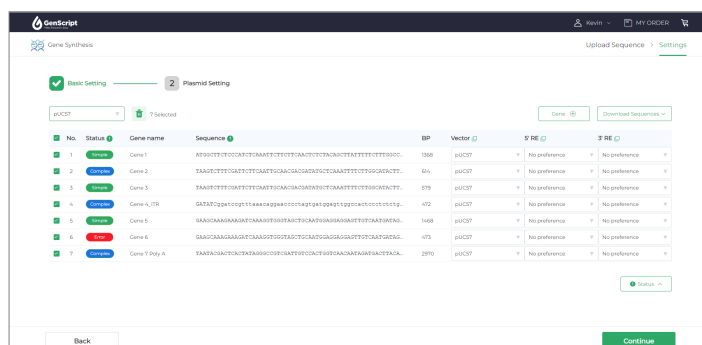


Figure 5: Sequence status. Once the vector and cloning method are determined, the sequences will be automatically analyzed for synthesis complexity and error-checking.



Figure 6: Sequence editing page. Click on the sequences to jump into sequence editing page.



Figure 7: Codon optimization page. Select the region, host organism, and restriction sites to keep and/or avoid prior to perform codon optimization.



Figure 8: Codon optimization result. Check the codon optimization result, you can continue the order with your original sequence or optimized sequence. You can also download the codon optimization report by clicking on the optimization report icon.

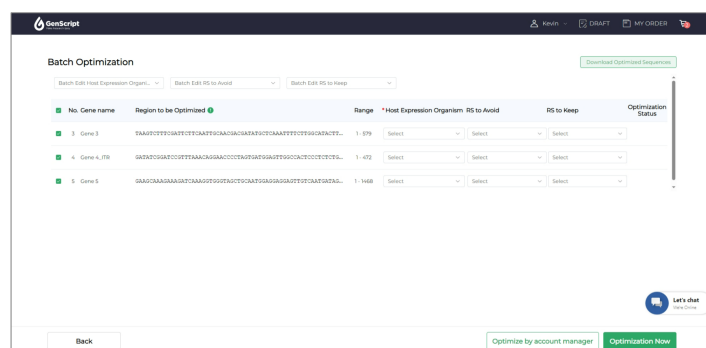
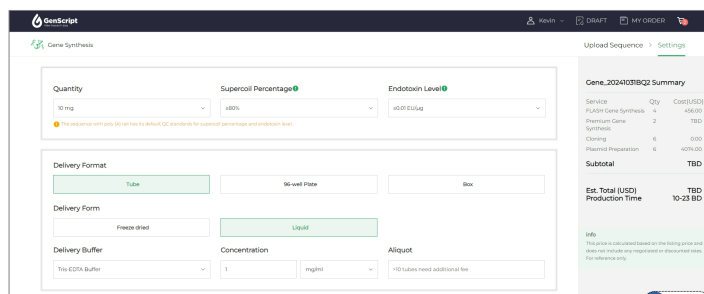


Figure 9: Batch codon optimization page. Batch select the region, host organism, and restriction sites to keep and/or avoid prior for batch codon optimization.

3. Plasmid Setup and Add-ons

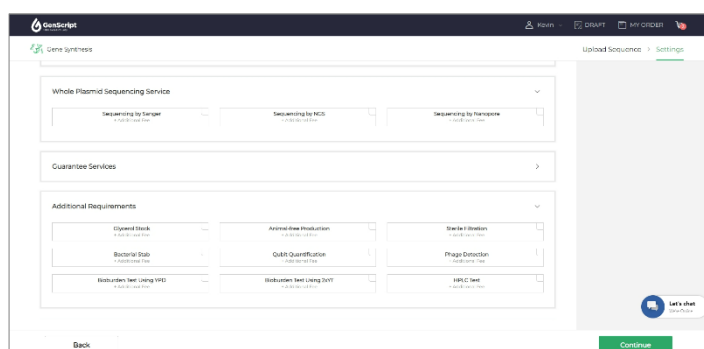
- The default plasmid DNA quantity is 4 micrograms for high-copy plasmids, but you can add buddled plasmid DNA preparation services to acquire high-quality gram-level plasmid DNA. Select the quantity, supercoil percentage and endotoxin level for your clonal genes based on your downstream application requirements (see **Figure 9**).
- Choose the delivery format and form for your plasmid DNA, and you can also customize the buffer and concentration (see **Figure 10**).
 ⚠ Please note that if your sequences contain Poly A structure or AAV ITR structure, Poly A guarantee and AAV ITR guarantee services are available to add to your order to ensure the complete and accurate Poly A and AAV ITR structure on your deliverables (see **Figure 11**).
- Add whole plasmid sequencing and other add-on services simply by clicking on those options (see **Figure 10**).
 ⚠ Please note that certain add-on services may require larger amount of plasmid DNA other than default 4 micrograms.



The interface shows settings for a plasmid DNA order. On the left, there are dropdown menus for Quantity (10 mg), Supercoil Percentage (x80%), and Endotoxin Level (x80% EndoLg). Below these are options for Delivery Format (Tube, 96-well Plate, Box) and Delivery Form (Freeze dried, Liquid). There are also fields for Delivery Buffer (Trio EDTA Buffer), Concentration (1 mg/ml), and Aliquot (100 µl). On the right, a 'Gene_2024020802 Summary' table lists services and costs.

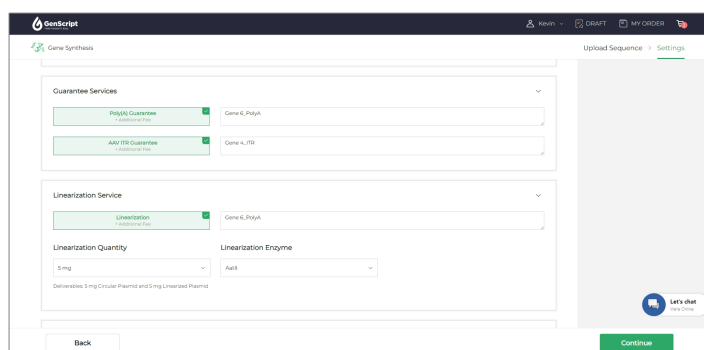
Service	Qty	Cost (USD)
PLASMID Gene Synthesis	4	450.00
Plasmid Gene Synthesis	2	TBD
Cloning	4	0.00
Plasmid Preparation	6	<0.01
Subtotal		TBD
Est. Total (USD)		TBD
Production Time		10-20 BD

Figure 9: Quantity, supercoil percentage, endotoxin level and delivery format & form settings. Depends on your application, choose the most suitable options of your plasmid DNA.



The interface shows the 'Whole Plasmid Sequencing Service' section with three dropdown menus: 'Sequencing by Target', 'Sequencing by T7E1', and 'Sequencing by Nanopore'. Below this is a 'Guarantee Services' section with a dropdown menu. The 'Additional Requirements' section includes checkboxes for 'Client Stock', 'Arrival date Production', 'Bacterial Strain', 'Quick Quarantine', 'Phage Detection', 'Bacterial test using VRE', 'Bacterial test using T4T', and 'EMC test'. A 'Continue' button is at the bottom right.

Figure 10: Whole plasmid sequencing and additional requirements. Add whole plasmid sequencing service and add-ons for your plasmid DNA.



The interface shows the 'Guarantee Services' section with checkboxes for 'Poly(A) Guarantee' and 'AAV ITR Guarantee'. Below this is the 'Linearization Service' section with a dropdown menu for 'Linearization' and a 'Linearization Quantity' dropdown menu. A 'Continue' button is at the bottom right.

Figure 11: Poly (A) guarantee and AAV ITR guarantee add-on services. Add guarantee services to ensure the complete and accurate structure in your sequences. You can also add plasmid linearization service for the sequences added poly (A) guarantee.

4. Order or Get a Quote

- Check the order summary, available coupons and pricing agreement contracts, turnaround time, and estimated cost that showing on the right side of the page (see **Figure 12**).
- Add new addresses or edit existing ones, by clicking "Management" next to the shipping address and billing address. For the shipping option, click "Edit" to select your preferred choice (see **Figure 12**).
- Once you have everything set up, click "Order Now" to place your order, or if you prefer to receive a detailed quote first, click "Get Quote" to request a quote for your order.
- You can use the existing payment method, or add new a new payment method on the payment page (see **Figure 13**).
- To keep track of your ongoing and past orders as well as quotations, click "My Order" at the top of the page (see **Figure 14**).

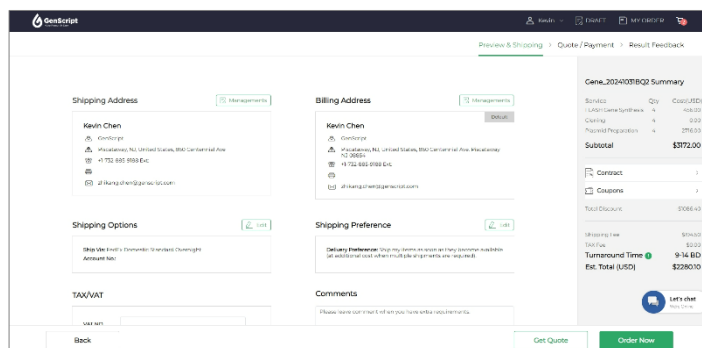


Figure 12 shows the GenScript checkout page. It includes sections for Shipping Address, Billing Address, Shipping Options, and Shipping Preference. A summary table on the right lists items and their costs.

Service	Qty	Cost (USD)
GenScript Synthesis	4	448.00
Cloning	4	8.00
Protein Purification	4	276.00
Subtotal		\$732.00
Total Discount		\$106.43
Shipping Fee		\$14.00
Sub Total		\$639.57
Turnaround Time		9-14 BD
Est. Total (USD)		\$2280.10

Figure 12: Available pricing agreement contracts and coupons, shipping address, billing address and shipping options. Select one available contract or coupon. Fill the shipping address and billing address, as well as select the shipping options for your order.

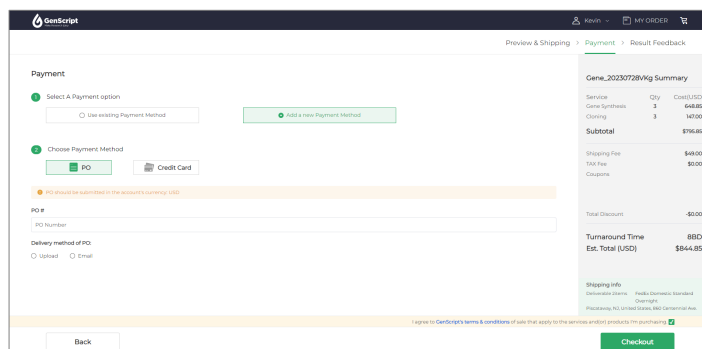


Figure 13 shows the GenScript payment page. It includes sections for Payment, Select A Payment option, and Choose Payment Method. A summary table on the right lists items and their costs.

Service	Qty	Cost (USD)
GenScript Synthesis	3	648.85
Cloning	3	14.00
Subtotal		\$762.85
Shipping Fee		\$40.00
Sub Total		\$802.85
Total Discount		\$0.00
Turnaround Time		8BD
Est. Total (USD)		\$844.85

Figure 13: Payment. Use existing payment method, or add new a new payment method.

GenScript

REAGENT SERVICESBIOINFORMATICSPRODUCTSAPPLICATIONSRESOURCESGROUP

Kevin Chen

Account Manager

Home

Projects

Draft

Quotation

Order

Review Request

FAQ

Vectors

Clones

Group

Account

Coupons & Rewards

Resources

Order

All Status

All Types of Orders

2023-07-26

2023-07-26

Order Name, ID or Gene, Plasmid, Cligle Name

View

Clear

Tip: If you need to find orders at a specific time, you can use the time filter.

Tip: If you have any questions about your orders please click the link and add the number, you'll get help from our customer service.

Click here for contact information.

Order ID	Type	Operation	Customized Name	PO	Order Date	Production Status	
GEN1782828	Gene		Order - customized name here		2023-07-26 9:10:23	Quote	
GEN1782829	Gene		Order - customized name here		2023-07-26 9:10:40	Quote	
GEN1782830	Gene		Order - customized name here		2023-07-26 9:10:57	Quote	
GEN1782831	Gene		Order - customized name here		2023-07-26 9:11:14	Quote	
GEN1782832	Gene		Order - customized name here		2023-07-26 9:11:31	Quote	
GEN1782833	Gene		Order - customized name here		2023-07-26 9:11:48	Quote	
GEN1782834	Gene		Order - customized name here		2023-07-26 9:12:05	Quote	
GEN1782835	Gene		Order - customized name here		2023-07-26 9:12:22	Quote	
GEN1782836	Gene		Order - customized name here		2023-07-26 9:12:39	Quote	
GEN1782837	Gene		Order - customized name here		2023-07-26 9:12:56	Quote	
GEN1782838	Gene		Order - customized name here		2023-07-26 9:13:13	Quote	
GEN1782839	Gene		Order - customized name here		2023-07-26 9:13:30	Quote	
GEN1782840	Gene		Order - customized name here		2023-07-26 9:13:47	Quote	
GEN1782841	Gene		Order - customized name here		2023-07-26 9:14:04	Quote	
GEN1782842	Gene		Order - customized name here		2023-07-26 9:14:21	Quote	
GEN1782843	Gene		Order - customized name here		2023-07-26 9:14:38	Quote	
GEN1782844	Gene		Order - customized name here		2023-07-26 9:14:55	Quote	
GEN1782845	Gene		Order - customized name here		2023-07-26 9:15:12	Quote	
GEN1782846	Gene		Order - customized name here		2023-07-26 9:15:29	Quote	
GEN1782847	Gene		Order - customized name here		2023-07-26 9:15:46	Quote	
GEN1782848	Gene		Order - customized name here		2023-07-26 9:16:03	Quote	
GEN1782849	Gene		Order - customized name here		2023-07-26 9:16:20	Quote	
GEN1782850	Gene		Order - customized name here		2023-07-26 9:16:37	Quote	
GEN1782851	Gene		Order - customized name here		2023-07-26 9:16:54	Quote	
GEN1782852	Gene		Order - customized name here		2023-07-26 9:17:11	Quote	
GEN1782853	Gene		Order - customized name here		2023-07-26 9:17:28	Quote	
GEN1782854	Gene		Order - customized name here		2023-07-26 9:17:45	Quote	
GEN1782855	Gene		Order - customized name here		2023-07-26 9:18:02	Quote	
GEN1782856	Gene		Order - customized name here		2023-07-26 9:18:19	Quote	
GEN1782857	Gene		Order - customized name here		2023-07-26 9:18:36	Quote	
GEN1782858	Gene		Order - customized name here		2023-07-26 9:18:53	Quote	
GEN1782859	Gene		Order - customized name here		2023-07-26 9:19:10	Quote	
GEN1782860	Gene		Order - customized name here		2023-07-26 9:19:27	Quote	
GEN1782861	Gene		Order - customized name here		2023-07-26 9:19:44	Quote	
GEN1782862	Gene		Order - customized name here		2023-07-26 9:20:01	Quote	
GEN1782863	Gene		Order - customized name here		2023-07-26 9:20:18	Quote	
GEN1782864	Gene		Order - customized name here		2023-07-26 9:20:35	Quote	
GEN1782865	Gene		Order - customized name here		2023-07-26 9:20:52	Quote	
GEN1782866	Gene		Order - customized name here		2023-07-26 9:21:09	Quote	
GEN1782867	Gene		Order - customized name here		2023-07-26 9:21:26	Quote	
GEN1782868	Gene		Order - customized name here		2023-07-26 9:21:43	Quote	
GEN1782869	Gene		Order - customized name here		2023-07-26 9:22:00	Quote	
GEN1782870	Gene		Order - customized name here		2023-07-26 9:22:17	Quote	
GEN1782871	Gene		Order - customized name here		2023-07-26 9:22:34	Quote	
GEN1782872	Gene		Order - customized name here		2023-07-26 9:22:51	Quote	
GEN1782873	Gene		Order - customized name here		2023-07-26 9:23:08	Quote	
GEN1782874	Gene		Order - customized name here		2023-07-26 9:23:25	Quote	
GEN1782875	Gene		Order - customized name here		2023-07-26 9:23:42	Quote	
GEN1782876	Gene		Order - customized name here		2023-07-26 9:23:59	Quote	
GEN1782877	Gene		Order - customized name here		2023-07-26 9:24:16	Quote	
GEN1782878	Gene		Order - customized name here		2023-07-26 9:24:33	Quote	
GEN1782879	Gene		Order - customized name here		2023-07-26 9:24:50	Quote	
GEN1782880	Gene		Order - customized name here		2023-07-26 9:25:07	Quote	
GEN1782881	Gene		Order - customized name here		2023-07-26 9:25:24	Quote	
GEN1782882	Gene		Order - customized name here		2023-07-26 9:25:41	Quote	
GEN1782883	Gene		Order - customized name here		2023-07-26 9:25:58	Quote	
GEN1782884	Gene		Order - customized name here		2023-07-26 9:26:15	Quote	
GEN1782885	Gene		Order - customized name here		2023-07-26 9:26:32	Quote	
GEN1782886	Gene		Order - customized name here		2023-07-26 9:26:49	Quote	
GEN1782887	Gene		Order - customized name here		2023-07-26 9:27:06	Quote	
GEN1782888	Gene		Order - customized name here		2023-07-26 9:27:23	Quote	
GEN1782889	Gene		Order - customized name here		2023-07-26 9:27:40	Quote	
GEN1782890	Gene		Order - customized name here		2023-07-26 9:27:57	Quote	
GEN1782891	Gene		Order - customized name here		2023-07-26 9:28:14	Quote	
GEN1782892	Gene		Order - customized name here		2023-07-26 9:28:31	Quote	
GEN1782893	Gene		Order - customized name here		2023-07-26 9:28:48	Quote	
GEN1782894	Gene		Order - customized name here		2023-07-26 9:29:05	Quote	
GEN1782895	Gene		Order - customized name here		2023-07-26 9:29:22	Quote	
GEN1782896	Gene		Order - customized name here		2023-07-26 9:29:39	Quote	
GEN1782897	Gene		Order - customized name here		2023-07-26 9:29:56	Quote	
GEN1782898	Gene		Order - customized name here		2023-07-26 9:30:13	Quote	
GEN1782899	Gene		Order - customized name here		2023-07-26 9:30:30	Quote	
GEN1782900	Gene		Order - customized name here		2023-07-26 9:30:47	Quote	
GEN1782901	Gene		Order - customized name here		2023-07-26 9:31:04	Quote	
GEN1782902	Gene		Order - customized name here		2023-07-26 9:31:21	Quote	
GEN1782903	Gene		Order - customized name here		2023-07-26 9:31:38	Quote	
GEN1782904	Gene		Order - customized name here		2023-07-26 9:31:55	Quote	
GEN1782905	Gene		Order - customized name here		2023-07-26 9:32:12	Quote	
GEN1782906	Gene		Order - customized name here		2023-07-26 9:32:29	Quote	
GEN1782907	Gene		Order - customized name here		2023-07-26 9:32:46	Quote	
GEN1782908	Gene		Order - customized name here		2023-07-26 9:33:03	Quote	
GEN1782909	Gene		Order - customized name here		2023-07-26 9:33:20	Quote	
GEN1782910	Gene		Order - customized name here		2023-07-26 9:33:37	Quote	
GEN1782911	Gene		Order - customized name here		2023-07-26 9:33:54	Quote	
GEN1782912	Gene		Order - customized name here		2023-07-26 9:34:11	Quote	
GEN1782913	Gene		Order - customized name here		2023-07-26 9:34:28	Quote	
GEN1782914	Gene		Order - customized name here		2023-07-26 9:34:45	Quote	
GEN1782915	Gene		Order - customized name here		2023-07-26 9:35:02	Quote	
GEN1782916	Gene		Order - customized name here		2023-07-26 9:35:19	Quote	
GEN1782917	Gene		Order - customized name here		2023-07-26 9:35:36	Quote	
GEN1782918	Gene		Order - customized name here		2023-07-26 9:35:53	Quote	
GEN1782919	Gene		Order - customized name here		2023-07-26 9:36:10	Quote	
GEN1782920	Gene		Order - customized name here		2023-07-26 9:36:27	Quote	
GEN1782921	Gene		Order - customized name here		2023-07-26 9:36:44	Quote	
GEN1782922	Gene		Order - customized name here		2023-07-26 9:37:01	Quote	
GEN1782923	Gene		Order - customized name here		2023-07-26 9:37:18	Quote	
GEN1782924	Gene		Order - customized name here		2023-07-26 9:37:35	Quote	
GEN1782925	Gene		Order - customized name here		2023-07-26 9:37:52	Quote	
GEN1782926	Gene		Order - customized name here		2023-07-26 9:38:09	Quote	
GEN1782927	Gene		Order - customized name here		2023-07-26 9:38:26	Quote	
GEN1782928	Gene		Order - customized name here		2023-07-26 9:38:43	Quote	
GEN1782929	Gene		Order - customized name here		2023-07-26 9:39:00	Quote	
GEN1782930	Gene		Order - customized name here		2023-07-26 9:39:17	Quote	
GEN1782931	Gene		Order - customized name here		2023-07-26 9:39:34	Quote	
GEN1782932	Gene		Order - customized name here		2023-07-26 9:39:51	Quote	
GEN1782933	Gene		Order - customized name here		2023-07-26 9:40:08	Quote	
GEN1782934	Gene		Order - customized name here		2023-07-26 9:40:25	Quote	
GEN1782935	Gene		Order - customized name here		2023-07-26 9:40:42	Quote	
GEN1782936	Gene		Order - customized name here		2023-07-26 9:40:59	Quote	
GEN1782937	Gene		Order - customized name here		2023-07-26 9:41:16	Quote	
GEN1782938	Gene		Order - customized name here		2023-07-26 9:41:33	Quote	
GEN1782939	Gene		Order - customized name here		2023-07-26 9:41:50	Quote	
GEN1782940	Gene		Order - customized name here		2023-07-26 9:42:07	Quote	
GEN1782941	Gene		Order - customized name here		2023-07-26 9:42:24	Quote	
GEN1782942	Gene		Order - customized name here		2023-07-26 9:42:41	Quote	
GEN1782943	Gene		Order - customized name here		2023-07-26 9:42:58	Quote	
GEN1782944	Gene		Order - customized name here		2023-07-26 9:43:15	Quote	
GEN1782945	Gene		Order - customized name here		2023-07-26 9:43:32	Quote	
GEN1782946	Gene		Order - customized name here		2023-07-26 9:43:49	Quote	
GEN1782947	Gene		Order - customized name here		2023-07-26 9:44:06	Quote	
GEN1782948	Gene		Order - customized name here		2023-07-26 9:44:23	Quote	
GEN1782949	Gene		Order - customized name here		2023-07-26 9:44:40	Quote	
GEN1782950	Gene		Order - customized name here		2023-07-26 9:44:57	Quote	
GEN1782951	Gene		Order - customized name here		2023-07-26 9:45:14	Quote	
GEN1782952	Gene		Order - customized name here		2023-07-26 9:45:31	Quote	
GEN1782953	Gene		Order - customized name here		2023-07-26 9:45:48	Quote	
GEN1782954	Gene		Order - customized name here		2023-07-26 9:46:05	Quote	
GEN1782955	Gene		Order - customized name here		2023-07-26 9:46:22	Quote	
GEN1782956	Gene		Order - customized name here		2023-07-26 9:46:39	Quote	
GEN1782957	Gene		Order - customized name here		2023-07-26 9:46:56	Quote	
GEN1782958	Gene		Order - customized name here		2023-07-26 9:47:13	Quote	
GEN1782959	Gene		Order - customized name here		2023-07-26 9:47:30	Quote	
GEN1782960	Gene		Order - customized name here		2023-07-26 9:47:47	Quote	
GEN1782961	Gene		Order - customized name here		2023-07-26 9:48:04	Quote	
GEN1782962	Gene		Order - customized name here		2023-07-26 9:48:21	Quote	
GEN1782963	Gene		Order - customized name here		2023-07-26 9:48:38	Quote	
GEN1782964	Gene		Order - customized name here		2023-07-26 9:48:55	Quote	
GEN1782965	Gene		Order - customized name here		2023-07-26 9:49:12	Quote	
GEN1782966	Gene		Order - customized name here		2023-07-26 9:49:29	Quote	
GEN1782967	Gene		Order - customized name here		2023-07-26 9:49:46	Quote	
GEN1782968	Gene		Order - customized name here		2023-07-26 9:50:03	Quote	
GEN1782969	Gene		Order - customized name here		2023-07-26 9:50:20	Quote	
GEN1782970	Gene		Order - customized name here		2023-07-26 9:50:37	Quote	
GEN1782971	Gene		Order - customized name here		2023-07-26 9:50:54	Quote	
GEN1782972	Gene		Order - customized name here		2023-07-26 9:51:11	Quote	
GEN1782973	Gene		Order - customized name here		2023-07-26 9:51:28	Quote	
GEN1782974	Gene		Order - customized name here		2023-07-26 9:51:45	Quote	
GEN1782975	Gene		Order - customized name here		2023-07-26 9:52:02	Quote	
GEN1782976	Gene		Order - customized name here		2023-07-26 9:52:19	Quote	
GEN1782977	Gene		Order - customized name here		2023-07-26 9:52:36	Quote	
GEN1782978	Gene		Order - customized name here		2023-07-26 9:52:53	Quote	
GEN1782979	Gene		Order - customized name here		2023-07-26 9:53:10	Quote	
GEN1782980	Gene		Order - customized name here		2023-07-26 9:53:27	Quote	
GEN1782981	Gene		Order - customized name here		2023-07-26 9:53:44	Quote	
GEN1782982	Gene		Order - customized name here		2023-07-26 9:54:01	Quote	
GEN1782983	Gene		Order - customized name here		2023-07-26 9:54:18	Quote	
GEN1782984	Gene		Order - customized name here		2023-07-26 9:54:35	Quote	
GEN1782985	Gene		Order - customized name here		2023-07-26 9:54:52	Quote	
GEN1782986	Gene		Order - customized name here		2023-07-26 9:55:09	Quote	
GEN1782987	Gene		Order - customized name here		2023-07-26 9:55:26	Quote	
GEN1782988	Gene		Order - customized name here		2023-07-26 9:55:43	Quote	
GEN1782989	Gene		Order - customized name here		2023-07-26 9:56:00	Quote	
GEN1782990	Gene	</					