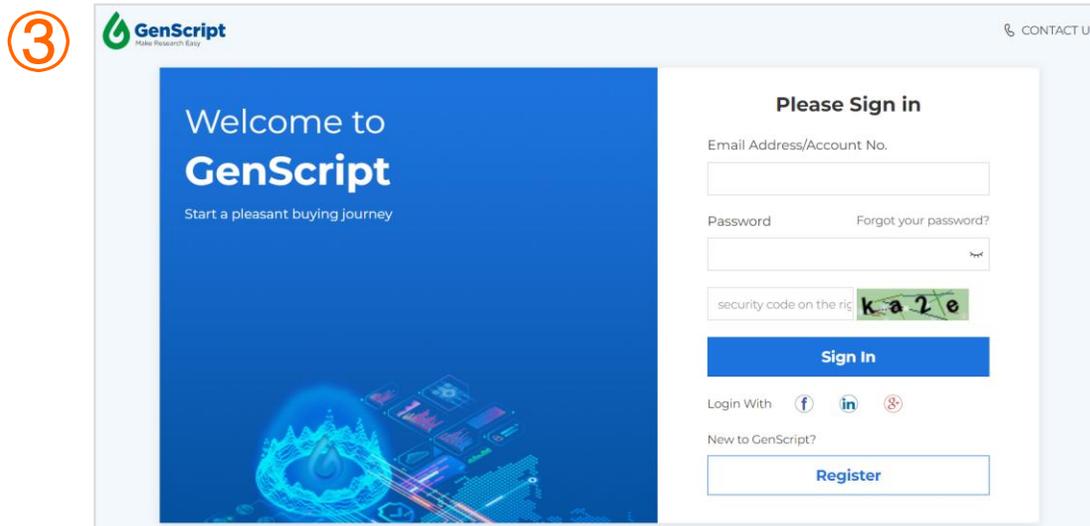
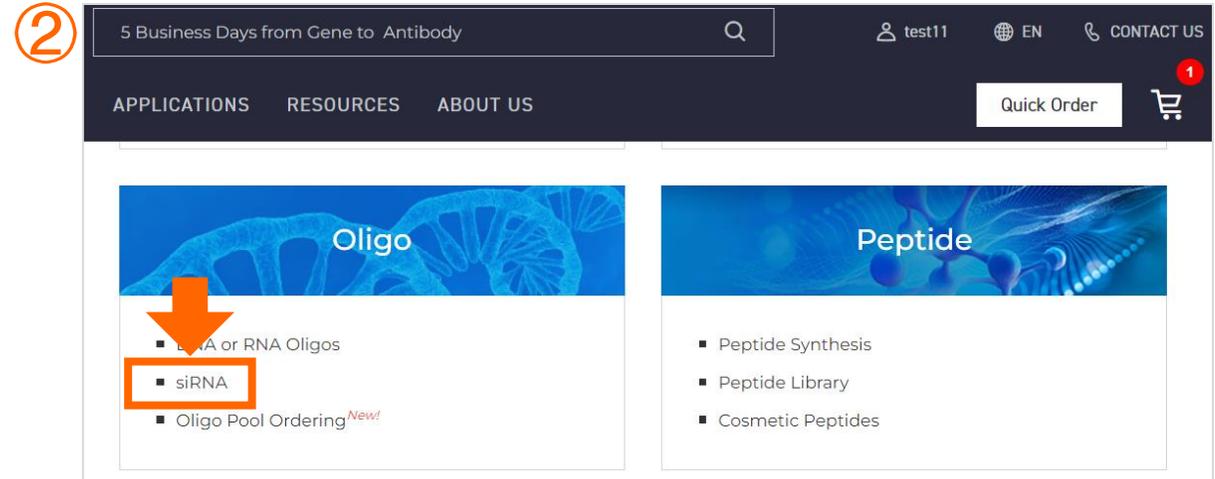
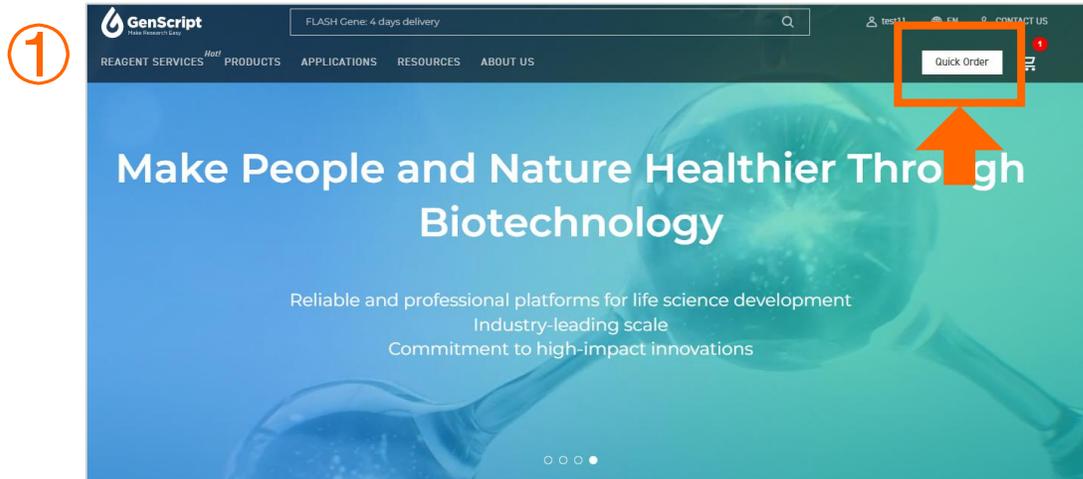


# User Guide for Custom siRNA Online Ordering

Please visit the [GenScript website](#). ① Select “Quick Order.” ② Select “siRNA” under Oligo services. ③ Register an account or log in.



# Step 1. Enter order requirements

① In **Order Information**: specify annealing preference and delivery format for all items in your order.

② In **Order Method: For Order by item**: Enter the desired quantity of siRNA.

**For bulk Order**: Download and complete the bulk order template, then upload it.

siRNA Ordering (\* Required Fields) 1 Information > 2 Cart > 3 Confirm Order > 4 Result Feedback

You have the latest draft to restore. Do you agree?  Auto save to draft: 02:45

## ① Order Information

Order Name:

\* Annealing:  ▼

\* Delivery Format:  ▼

\* Container:  ▼

96-well plate display area  
(shown when 96-well plates are selected)

## ② Order Method (choose 1 of 2)

**Order by item:** Number of siRNA:

**Bulk Order:**

## Step 2. Enter siRNA sequences and specifications

- ① Enter sense and antisense sequences manually *or* copy/paste them directly from Excel into the ordering form.
  - Use 'mA,' 'mU,' 'mC,' and 'mG' for 2'-O-methylation bases.
  - Add '\*' between two bases for phosphorothioate modifications.
  - Use 'dT' for T (Thymidine).
- ② Select additional specifications like purification level and quantity for each items.
- ③ Use the functional buttons for all sequences as needed.
- ④ Select positive and negative controls in the add-on service if needed.

③

Automatically fill in antisense strand based on sense strand Add full-length 2'-O-Methylation modification Add full-length phosphorothioate modification T to U Clear Table

Form instructions: Bulk copy/paste from an Excel sheet here ▼ Number in sequence ☑ Fill according to first row ☑ Clear all values

	*siRNA Name ↓ ☑	*Sense Sequence 📄 ☑ ?	*Anti-sense Sequence 📄 ☑ ?	5' Modification 📄 ☑	3' Modification 📄 ☑	Internal Modification ☑	*Purification 📄 ☑	*Unit 📄 ☑	*Quantity 📄 ☑	*Aliquoting Into 📄 ☑	Price	Opti
1				+	+		+ RNase free HPLC	nmol				👁️ 🗑️
2				+	+		+ RNase free HPLC	nmol				👁️ 🗑️
3				+	+		+ RNase free HPLC	nmol				👁️ 🗑️
4				+	+		+ RNase free HPLC	nmol				👁️ 🗑️
5				+	+		+ RNase free HPLC	nmol				👁️ 🗑️

Add  rows Apply  Confidential sequence ?

If you have any special requirements, for example adding any modifications to the RNA sequence, please describe them here.

④

Add-on Items (positive and negative controls)

# Step 3. Add modifications

## 5'/3' Modifications

- 1 Click a "+" icon in the 5' or 3' Modification columns.
- 2 Choose sense or antisense sequence to modify.
- 3 Use the search function to quickly find your modification.
- 4 Choose your modification and click "Confirm."

The screenshot shows the '5'/3' Modification' interface. A table with columns for 'Sense Sequence', 'Anti-sense Sequence', and '5' Modification' is visible. A '+' icon in the '5' Modification' column is circled with a '1'. A dropdown menu for 'Modification Position' is set to 'Sense Sequence' and circled with a '2'. A search box for 'Terminal Modification Name' is circled with a '3'. A list of modifications is shown, with '5' Phosphate' circled with a '4'. A 'Confirm' button is at the bottom right.

## Internal Modifications

- 1 Click a "+" icon in the Internal Modification column.
- 2 Choose sense or antisense sequence to modify.
- 3 Use the search function to quickly find your modification.
- 4 Place the cursor where you want the modification.
- 5 Select modification from the list and click "Add."

The screenshot shows the 'Internal Modification' interface. A dropdown menu for 'Modification Position' is set to 'Sense Sequence' and circled with a '2'. A search box for 'Internal Modification Name' is circled with a '3'. A text input field for 'Sense Sequence' is circled with a '4'. A list of modifications is shown, with the 'Action' column containing 'Add' buttons circled with a '5'. A '+' icon in the 'Internal Modification' column of a table is circled with a '1'. A 'Confirm' button is at the bottom right.

# Step 4. Confirm price and production time

- ① Verify the list price for each item.
- ② Select a contract code or coupon if applicable.
- ③ Review the estimated production time and order total.
- ④ Click “Add to Cart” and proceed to checkout.

Add full-length 2'-O-Methylation modification

Add full-length phosphorothioate modification

T to U

Clear Table

Guidance

Form instructions: Bulk copy/paste from an Excel sheet here ↓ Number in sequence 📄 Fill according to first row 🗑️ Clear all values

	* siRNA Name ↓ 📄	* Sense Sequence 📄 🗑️ ⓘ	* Anti-sense Sequence 📄 🗑️ ⓘ	5' Modification 📄 🗑️	3' Modification 📄 🗑️	Internal Modification 🗑️	* Purification 📄 🗑️	* Unit 📄 🗑️	* Quantity 📄 🗑️	* Aliquoting Into 📄 🗑️	Price	Opti
1	1	aucgaucgaucgaucgaucgaucg	gcaugcaugcaugcaugcaugcau	+	+	+	RNase free HPLC ▾	nmol	2	1	\$ 59.00	🗑️
2	2	gcaugcaugcaugcaugcaugcau	aucgaucgaucgaucgaucgaucg	+	+	+	RNase free HPLC ▾	nmol	10	1	\$ 123.00	🗑️
3				+	+	+	RNase free HPLC ▾	nmol	+			🗑️
4				+	+	+	RNase free HPLC ▾	nmol	+			🗑️
5				+	+	+	RNase free HPLC ▾	nmol	+			🗑️

Add  rows Apply

Confidential sequence ⓘ

If you have any special requirements, for example adding any modifications to the RNA sequence, please describe them here.

③

②

Contract Save: \$0.00   
  Coupon Save: \$45.50  
▼ Discount Details

③

Estimated Time: 5 Business Days

Total Amount: \$182.00

Save Draft

④

Add To Cart