

Custom Small RNA Blocker – Request Form (Fillable)

Complete form and send it by e-mail to ngs@revvity.com. All fields required. Research Use Only

1) Customer & Order Details

Company name	<input type="text"/>		
Contact name	<input type="text"/>	Title	<input type="text"/>
Email	<input type="text"/>	Phone	<input type="text"/>

2) Project Overview

Goal:	Deplete specific miRNAs	Reduce tRNA or YRNA fragments
Other:	<input type="text"/>	

3) Biological Context

Species	<input type="text"/>				
Sample types (check all that apply):	Plasma	Serum	Tissue	Cell line	Plant
EVs	Other:	<input type="text"/>			

4) Targets To Block

How many targets to block (count) - 10 max	<input type="text"/>
--	----------------------

Design Note:

For each target, provide exact RNA sequence (5'→3') in FASTA format.

Example FASTA entry:

```
>mir-XYZ_hsa
AUGC UUACGAGUGACUUGA
```

Minimum Target Length: 20 nt
Maximum Target Length: 60 nt

FASTA entries (paste one per target):