

# ScreenToLead™ Oligos

## Supporting Candidate Screening

We have leveraged our expertise to design oligo packages with **defined parameters**, delivering the most suitable antisense oligonucleotide (ASO) sequences to support your **candidate screening** process.

### ScreenToLead™ Oligo Specifications

SPECIFICATION	DETAILS
Length	15 to 22 mers
Quantity	≥10 nmoles/oligo
Chemistries and Modifications	LNA®, 2'O-Me, 2'O-MOE, 2'-Fluoro RNA Full Phosphorothioate backbone
Format	Desalted, dried in single tubes
Quality Controls	Mass spectrometry + single quantification
Documentation	Technical data sheet
Optional Purification	HPLC
Min ordering quantity	20 oligos per set
Average delivery	2 – 3 weeks

### ScreenToLead™ Oligo Packages

CAT #	CHEMISTRY	SPECIFICATIONS	PRICE			
SC-STL-Full	Up to <b>Full modified</b>	ScreenToLead™ Oligos - Up to a fully modified 15-22-mers antisense oligonucleotide	179 EUR	179 CHF	160 GBP	202 USD
SC-STL-GAP	<b>Gapmer</b>	ScreenToLead™ Oligos - 15-22-mers Gapmer antisense oligonucleotide - Maximum 10 modifications	105 EUR	105 CHF	94 GBP	119 USD

### Get Started easily

Send your sequences to [oligo.centre@eurogentec.com](mailto:oligo.centre@eurogentec.com) and receive a quote.

## State-of-the-art equipment

Our laboratories, equipped with state-of-the-art **Shasta synthesizers**, enable the **parallel synthesis of 5 to 192 oligonucleotides**, with the capacity to support more **extensive projects** upon request.

## Fast turnaround time production

We produce custom ASO sequences with a **fast turnaround time** to **accelerate** our customers' research and development **timelines** without compromising quality.

## Continued support beyond early stages

Our support and manufacturing capabilities extend beyond discovery and screening, offering a **continuum of services** from research to clinical. Customers can rely on a **single, trusted partner** throughout their **therapeutic development** journey.



With **40+** years of experience, we also provide **tailored solutions** for custom **antisense** oligonucleotide (ASO).



**40 years**  
of experience



**Hundreds**  
of sequences/  
order



Variety of  
**Modifications**



**Continuum**  
up to clinical

SCAN THIS CODE AND

**Discover our ASO offer**

