



FREQUENTLY ASKED QUESTIONS

Sigma Genosys and Sigma Proligo become ONE!

Sigma Genosys and Sigma Proligo become ONE!! Sigma-Aldrich will be launching a new online order system for Sigma Life Science custom oligonucleotides and peptides during September 2008. At the ONE website you will be able to order all your familiar custom products and much, much, much more...

1. *Will my custom oligonucleotide product selection change?* **Yes. For both our Sigma Genosys and Sigma Proligo customers there will be a range of new product selections.**

Standard Oligos

- Standard oligos (DNA and RNA) are ordered by manufacturing scale (not OD).
- Each manufacturing scale has a minimum yield guarantee based on purification, length and modifications. Please see Table 1 further information on ordering scales and guaranteed yield.
- All manufacturing scales, except 0.025 μ mol are full synthesis yield and have a minimum guaranteed yield as indicated in Table 1.
- For the 0.025 μ mol scale the yield is fixed at 3OD.
- Guaranteed Oligos™, Mass Oligos™ and Premium Oligos™ will be discontinued.
- E@sy Oligos™ (3OD, 100 μ M, Fixed price for 15-30 bases only) will be available as a standard product offering.
- 5', 3' and internal modifications are ordered by manufacturing scale.

Probes, Molecular Beacons and siRNA oligos

- Dual Labeled Fluorescent Probes (DLFPs), Lightcycler™ Probes, Scorpion™ Probes, Molecular Beacons and siRNA will be ordered by OD.
- DLFP Probes and Molecular Beacons will be available in the following scales - 1OD, 3OD, 5OD, 10OD, 50OD and 100OD.

2. *What is the new format for standard RNA oligos?*
 - RNA oligos will be sold by manufacturing scale.
3. *Can I order my standard oligos in solution?* **Yes.**
 - We can provide your oligos in solution (concentration adjusted).
4. *Since I do not place my oligo orders online, will I be impacted by the product alignment changes?* **Yes.**
 - Changes to the product portfolio will impact all Sigma Proligo and Sigma Genosys customers, even if you do not order online. All email orders must be submitted in the new format. Templates for the new format will be available online via the new online ordering website.

5. *Will pricing change? Yes.*

- Due to the consolidation of the product ranges, almost all pricing will have changed, however for most standard oligos (less than 50mer, desalted, no modifications), there should be minimal change to your current price. Price changes primarily affect modifications, specialty primers and purified oligos.
- There is now no additional charges for long oligos or wobbles.
- Pricing for standard DNA and RNA oligos is determined by the following formula:
$$(\text{number of bases} \times \text{price per base}) + \text{Purification Charge} + \text{Modification Charges (if requested)} = \text{Price per oligo}$$
- Please see Table 1 below for pricing on our most common oligos.

6. *Is there a change to the amount I pay for freight? Yes.*

- As from the 1st September 2008, if you order online you will receive a 20% discount off our standard freight fee of \$15 for Australia and \$20 for New Zealand.
- Freight for online orders will be \$12 for Australia and \$16 for New Zealand.

7. *Can I order items from the Sigma-Aldrich range of catalogue products at the same time as I order my oligos? Yes.*

- You can now order both custom oligos and peptides and other Sigma-Aldrich catalogue products on the one order. Please note, orders are not consolidated and will be shipped separately. Separate shipping assures that your order is not delayed. As a consequence, separate delivery charges will apply to both components (custom products and catalogue products) of the order.

8. *When do the changes take affect?*

- All product format and pricing changes will take affect on August 25th 2008.
- Pricing and format changes will be implemented for all our customers, regardless of how you order your oligos ie either via the existing Sigma Genosys or Sigma Proligo website, via the new website or email.
- For the most accurate information on product specifications and pricing, we recommend that you switch to our new site as soon as you are sent your login and password details. Look out for your email during the first two weeks of September.

Table 1 – Comparison of existing Sigma Genosys and Sigma Proligo product formats and the new Sigma Life Science product format

| Existing Product Format | | | | New Product Format | | |
|-----------------------------|-----------------------------|--------|---------------------------------|---------------------------------|------------|-----------------------|
| | Scale | Yield | Price (AUD) / Base | Scale | Yield | Price (AUD) / Base |
| Sigma Proligo | Gauranteed Oligos | | | DNA Oligos | | |
| | 2OD | 2OD | \$0.90 | 0.025µmol | 3OD | \$0.90 |
| | 5OD | 5OD | \$0.95 | 0.05µmol | > 5OD | \$0.95 |
| | 10OD | 10OD | \$1.85 | 0.20µmol | >12OD | \$2.45 |
| | 50OD | 50OD | \$5.50 | 1µmol | >40OD | \$4.45 |
| | E@sy Oligos | | | E@sy Oligos | | |
| | 3OD | 3OD | \$23/oligo | 3OD | 3OD | \$23/oligo |
| | Premium Oligos (RPC) | | | Base Price + RP1 Charge | | |
| | 2OD | 2OD | \$1.35 | 0.025µmol | NA | NA |
| | 4OD | 4OD | \$2.35 | 0.05µmol | > 1OD | \$15 |
| | NA | NA | NA | 0.20µmol | > 3OD | \$15 |
| | NA | NA | NA | 1µmol | > 12OD | \$25 |
| | PAGE Purified Oligos | | | Base Price + PAGE Charge | | |
| | NA | NA | NA | 0.05µmol | > 0.5OD | \$75 |
| | 1OD | 1OD | \$3.40 | 0.20µmol | > 1OD | \$75 |
| NA | NA | NA | 1µmol | > 5OD | \$88 | |
| HPLC Purified Oligos | | | Base Price + HPLC Charge | | | |
| 2OD | 2OD | \$2.50 | 0.05µmol | > 1OD | \$75 | |
| 5OD | 5OD | \$4.60 | 0.20µmol | > 2.5OD | \$75 | |
| 10OD | 10OD | \$7.65 | 1µmol | >13OD | \$88 | |
| Sigma Genosys | NA | NA | NA | 0.025µmol | 3OD | \$0.90 |
| | 0.05µmol | > 5OD | \$0.95 | 0.05µmol | > 5OD | \$0.95 |
| | 0.20µmol | >12OD | \$2.45 | 0.20µmol | >12OD | \$2.45 |
| | 1µmol | >40OD | \$4.45 | 1µmol | >40OD | \$4.45 |
| | Easy Oligos | | | Easy Oligo | | |
| NA | NA | NA | 3OD | 3OD | \$23/oligo | |

Note: Modification charges remain unchanged for Sigma Genosys customers

Note: Standard Oligos refer to oligos <50 bases, no modifications or wobbles