

MGB-Eclipse™ Probes

For any applications

License-free use

Patents on MGB Eclipse™ Probes have now expired. By buying your MGB Probes at Eurogentec, you are no longer limited by any licenses. You can now use our MGB-Probes for any applications in research, agro-food, vet and human management.

Custom Synthesis of MGB-Eclipse™ Probes

- From **8 to 30 bases**
- **6 - 20 - 50 nmol** delivered quantities
- **RP-HPLC** Purified
- **MALDI-TOF MS/UHPLC** quality controlled
- Delivered in **solution** (H₂O or TE); **Dried** or in **Dried Aliquotes** of 10 nmol

MGB Probes with Fluorescent Dyes covering all qPCR channels.



The Advantages of MGB probe technology



Specific

MGB Probes increase the T_m of a probe and offer a very high specificity toward its complementary sequence.



Sensitive

MGB Probes ensure the detection of low copy targets.



Efficient

MGB Probes are perfectly suited for patient management and the diagnostic of pathogens.

How to order on eurogentec.com?

Our MGB Configurator is the easiest way to order your MGB Probes.

1 Configure your MGB Probe on line

Define all the parameters for your MGB Probes: Sequence, Fluorescent Dyes and Quantity. You can also use the batch order file to order multiple MGB Probes at a time.

2 Benefit from 2 free primers

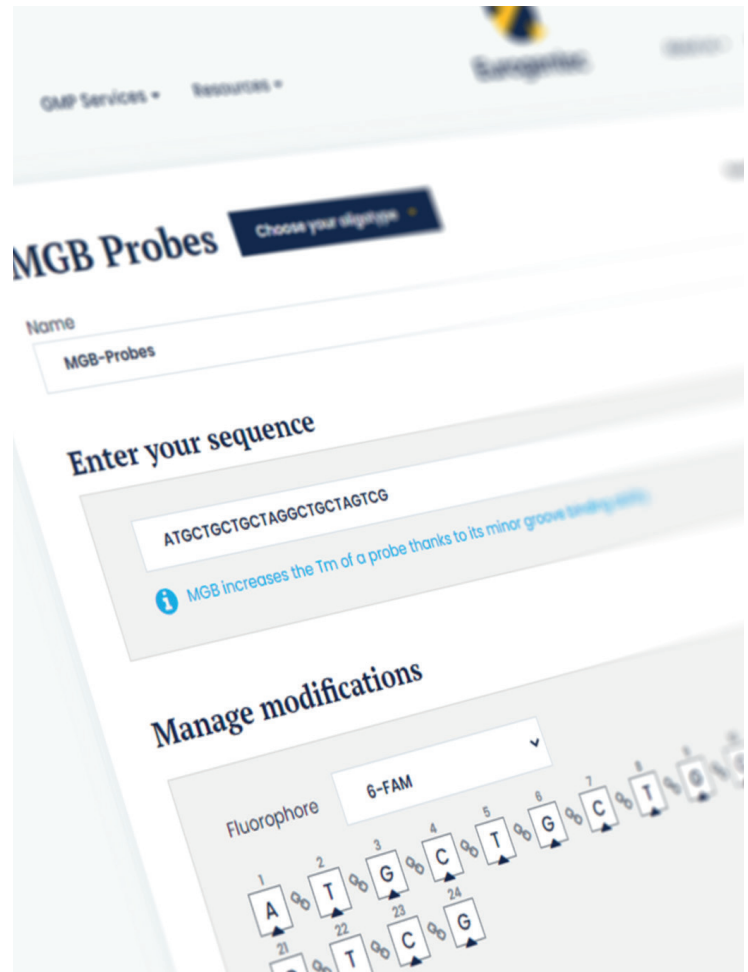
On the same screen, you can select up to 2 compatible free primers among your oligo list or configure new ones.

3 Add to Cart

Click on the Add to Cart button at the bottom of the MGB Configurator. Your probe and your free primer(s) will be added together in your cart and saved in your oligo list.

4 Confirm your order

Finalize your order



Why Choose MGB Probes from Eurogentec?

Expertise

We have been producing oligonucleotides for research and IVD applications **since 1985** and have expanded our portfolio with **ELITech Group's** MGB Probes in 2015.

Quality

We manufacture oligonucleotides from high quality raw materials. Our processes are **ISO 9001** and **ISO 13485** certified according to the oligo grade.

Partnership

We assist our customers in their **strategy** to commercialize their own diagnostic kits and offer them the **best approach** in compliance with the regulations.

Order on www.eurogentec.com