

BHQ[®] dual labeled qPCR Probes

for all applications

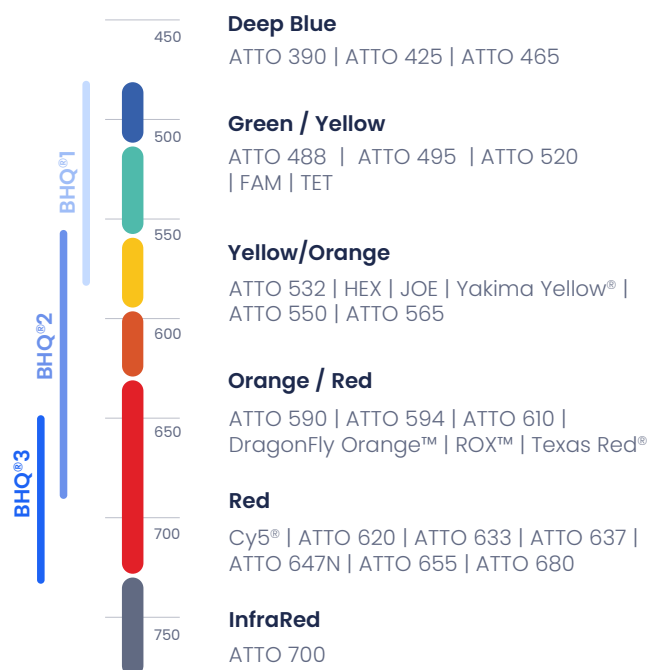
License free BHQ[®] from Eurogentec

We produce modified oligonucleotides and qPCR Probes labeled with **BHQ[®]-1**, **BHQ[®]-2** and **BHQ[®]-3**. Buying BHQ[®] labeled oligonucleotides from Eurogentec is **no longer limited by any licenses** for research, diagnostic and commercialization purposes.

Compatible with most of the popular reporter dyes

Choosing the right quencher for a qPCR probe is a critical aspect when developing an assay. To be efficient, the quencher absorption must cover the emission wavelength of the fluorescent dye. BHQs[®] have the advantage of **covering the full visible spectrum** and are compatible with most of the fluorescent dyes.

BHQ[®] Absorption Wavelength Range (in nm) and popular fluorescent dyes



Why use our BHQ labeled oligo?



License free for any Applications

We produce Custom labeled oligos with the original BHQ[®] quenchers **for any applications without license restriction.**



Broad absorption Range

BHQ[®] are compatible with a large range of fluorescent dyes as they absorb fluorescent emission **from 480 nm to 730 nm.**



Grade continuum from Research to Diagnostics

We produce Custom oligonucleotides in **various grades** according to your applications.

At your side from research to commercialization

We have a strong background in oligonucleotide production and offer suitable oligo grades from research to commercialization. Our scientific experts work in state-of-the-art facilities to deliver quality products in accordance with the customer's request.

Each oligo grade is accompanied with the appropriate documentation according to its usage.

Oligonucleotide Grades	Research	Track™	Pre-Diagnostic	Diagnostic
Process				
Dedicated Account Contact Person	Option	✓	✓	✓
Customized Fill & Finish	Option	Option	✓	✓
Quality Management				
ISO 9001 Certified Process	✓	✓	✓	✓
ISO 13485 Certified Process	-	-	✓	✓
Qualification/Validation [Equipment & Method]	-	Partial	✓	✓
Control				
Quantification	Single	Dual	Triple	Triple
Stringent QC Tests [validated]	-	✓	✓	✓
Traceability	Partial	Documented	Documented	Full Documented
Batch Record [Archived for 5 years]	-	-	Partial	Full
Classified Cleanrooms	-	-	✓	✓
Certificate of Analysis [CoA]	-	✓	✓	✓

BHQ® is a licensed trademark of Biosearch Technologies, Inc.

Order on www.eurogentec.com