

Access™ Dyes & Quenchers

From R&D to commercial use

Our **Access™ Dyes & Quenchers** consist of **IP-friendly** efficient reporters & quenchers for dual-labeled **qPCR and dPCR probes**.

The selection of dyes for **diagnostic tests** covers the **full visible spectrum**, from blue to far-red-shifted wavelengths.

BHQ® quenchers

Label your custom oligonucleotides and qPCR probes with the **original BHQ®** quencher dyes. It is now **license-free** at Eurogentec for research, diagnostic and commercial use. BHQ® quenchers cover **all qPCR channels** and are compatible with most of the fluorescent dyes.

MGB Eclipse® probes

Increase the specificity, sensitivity, and efficiency of your assay with our custom **license-free MGB Eclipse®** probes.

ATTO Dyes

Extend your possibilities by using robust and sensitive dyes covering the **full spectrum**. Moreover, ensure a steady supply by using **original molecules** for your commercial purposes.

A customizable and attractive offer



Friendly license

Friendly license plans for use in commercial applications



Broad range

Broad range of dyes including popular reporter/quencher pairs as well as proprietary molecules



Large synthesis scale

Small to **large synthesis scale** of dual-labeled custom oligonucleotides



Unique Partner

A **Unique Partner** for all your Nucleic Acid-Based Applications, with research, Track™ and diagnostic grade oligos

Find the most efficient dye and quencher pairs for qPCR Probes

Our selection of dyes for diagnostic tests covers all the full visible spectrum.

Channels	Dye and Quencher pairs	Dye Em max (nm)	Quencher Abs max (nm)
● Deep Blue	ATTO 390 / BHQ [®] -1	476	534
	ATTO 425 / BHQ [®] -1	485	534
● Green Yellow	FAM / BHQ [®] -1	520	534
	FAM / MGB-EDQ	520	342/545
● Yellow Orange	YY / BHQ [®] -1	549	534
	HEX / BHQ [®] -1	556	534
	HEX / MGB-EDQ	556	342/545
	ATTO 550 / BHQ [®] -2	576	579
● Orange Red	ROX / BHQ [®] -2	602	579
	ROX / MGB-EDQ	602	342/545
	ATTO 590 / BHQ [®] -2	622	579
● Red	ATTO 647 N / BHQ [®] -2	664	579
	ATTO 647 N / BHQ [®] -3	664	680
	HL647 / QXL [®] 670	675	668
● Infrared	HL750 / IRQXL [®]	778	772

Our fluorescent probes and quenchers are just **as effective in digital PCR.**

Configure your custom probes with our Access[™] Dyes & Quenchers

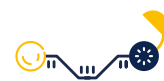


Double-Dye Probes

We provide classic and locked nucleic acid probes.

Specifications

Length: from 8 to 45 bases
Purification: HPLC
QC: MALDI-TOF MS + UPLC
Format: dried



MGB Probes

MGB increases the T_m of a probe thanks to its minor groove binding ability.

Specifications

Length: from 8 to 30 bases
Purification: HPLC
QC: MALDI-TOF MS + UHPLC
Format: In solution (H₂O or TE); dried or dried aliquoted in 10 nmol on request

Order on www.eurogentec.com