

ATTO Dyes for Probe Assays

For your research and commercial projects



15+ Licence free ATTO dyes

Discover the whole range on the back!

Ideal for commercial applications

Custom probes, labeled with IP-free ATTO Dyes from Eurogentec, are **ideal for commercial applications** such as human diagnostics, veterinary testing, and agro-food analysis. By choosing an ATTO Dye from the early discovery stage of your project, you ensure a **smooth path to commercialization without the burden of licensing**.

From Research to GMP-IVD grade

With extensive experience in oligonucleotide production, we provide oligo grades from research to in vitro diagnostics, each with the **required documentation according to its intended use**. Our team of scientific specialists operates in **cutting-edge facilities**, ensuring **high-quality products** tailored to meet each customer's specifications.

Why choose an ATTO dye for your probes?



Bright

ATTO dyes provide bright signals, enhancing the sensitivity and precision of your probes to ensure accuracy in nucleic acid-based assays.



Stable

ATTO dyes are photo- & thermo-stable, retaining their fluorescence intensity even at high temperatures, ensuring consistent and accurate results.



Steady supply

Our ATTO dyes are sourced directly from the original manufacturer, making them the reliable and secure choice for all your commercial projects.



Covering the full spectrum

Our selection of ATTO dyes ensures optimal performance across the full visible spectrum including infrared.

ATTO dyes offer a high-quality alternative to commonly used fluorophores. They can be combined with our IP-free quenchers to generate **double-dye probes** for commercial uses with **no IP**.

Our ATTO dyes are available for Double-Dye probes and MGB probes.

Channels	Dye	Dye Em max (nm)	Alternative to	IP-free Quenchers
Deep Blue	ATTO 390	476		BHQ [®] -1
	ATTO 425	485		BHQ [®] -1
Green Yellow	ATTO 465	509		BHQ [®] -1
	ATTO 488	520	FAM, Alexa Fluor 488	BHQ [®] -1, MGB-EDQ
	ATTO 520	538		BHQ [®] -1
Yellow Orange	ATTO 532	552	VIC [®] , HEX, Yakima Yellow [®]	BHQ [®] -1, MGB-EDQ
	ATTO 542	562		BHQ [®] -2
	ATTO 550	576	Cy3, Alexa Fluor 555	BHQ [®] -2
	ATTO 565	590	ROX, Cy3.5	BHQ [®] -2, MGB-EDQ
	ATTO RHO12	600		BHQ [®] -2
Orange Red	ATTO 590	622	Texas Red [®] , Alexa Fluor 594	BHQ [®] -2
	ATTO 594	626		BHQ [®] -2
Red	ATTO 633	651	LIZ [®]	BHQ [®] -2/-3
	ATTO 643	665		BHQ [®] -2/-3
	ATTO 647N	664	Cy5, Alexa Fluor 647	BHQ [®] -2/-3
	ATTO 655	680		BHQ [®] -2/-3
	ATTO 680	698	Cy5.5	BHQ [®] -2/-3
Infrared	ATTO 700	716	Cy5.5	BHQ [®] -3
	ATTO 740	763		IRQXL [®]

ATTO dyes are an ATTO-TEC technology



Order on www.eurogentec.com

