

# Spectral Properties and Quenching Range of Black Hole Quencher BHQ-1™, BHQ-2™ & BHQ-10™

Black Hole Quenchers™ (BHQs™) are the gold standard among quenchers in molecular biology, especially for quantitative PCR (qPCR) and other fluorescence-based assays. Unlike traditional quenchers, BHQs™ are true dark quenchers without native emission, which reduces background noise and increases detection sensitivity.

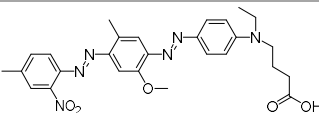
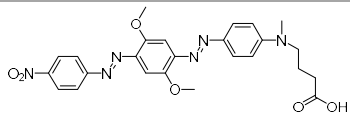
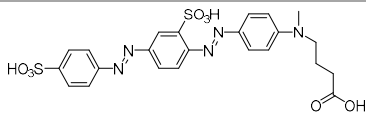
Jena Bioscience now offers BHQ-1™-, BHQ-2™- and BHQ-10™- labeled nucleotides to quench fluorescence across different spectral ranges:

BHQ-1 has an absorption maximum at 534 nm and a quenching range between 480 - 580 nm, making it suitable for quenching commonly used green and yellow fluorophores like Cy3, Fluorescein and Rhodamine. BHQ-2, with an absorption maximum at 579 nm, is used to quench in the range 560 – 670 nm and is most effective in quenching orange and red dyes such as Cy5, AF594 and Texas Red.

BHQ-10 is a water-soluble quencher with an absorption maximum at 615 nm and a quenching range of 480-550 nm. This makes BHQ-10 suitable for quenching blue and green fluorophores like AF488, Fluorescein and Rhodamine (Table 1).

**Table 1: Spectral properties and quenching range of Black Hole Quencher BHQ-1™, BHQ-2™ & BHQ-10™.**

**Please note:** It is important that the emission peak of the fluorescent dye and the absorption peak of the quencher have a large overlap range that ensures efficient resonant energy transfer and thus efficient quenching.

Dye	Quencher		
	BHQ-1™	BHQ-2™	BHQ-10™
	$\lambda_{exc}$ 534 nm $\epsilon$ 34.0 L mmol <sup>-1</sup> cm <sup>-1</sup>	$\lambda_{exc}$ 579 nm $\epsilon$ 38.0 L mmol <sup>-1</sup> cm <sup>-1</sup>	$\lambda_{exc}$ 516 nm $\epsilon$ 28.7 L mmol <sup>-1</sup> cm <sup>-1</sup>
			
<b>IR750</b>	-	-	-
<b>ATTO 680</b>	-	-	-
<b>IR680LT</b>	-	-	-
<b>ATTO 655</b>	-	-	-
<b>AF647</b>	-	-	-
<b>Cy5</b>	-	<b>x</b>	-
<b>ATTO 647N</b>	-	<b>x</b>	-
<b>ATTO 643</b>	-	<b>x</b>	-
<b>STAR RED</b>	-	<b>x</b>	-

# Spectral Properties and Quenching Range of Black Hole Quencher BHQ-1™, BHQ-2™ & BHQ-10™

<b>ATTO 594</b>	-	X	-
<b>AF594</b>	-	X	-
<b>Texas Red</b>	-	X	-
<b>STAR 580</b>	-	X	-
<b>AF555</b>	X	X	-
<b>ATTO 550</b>	X	X	-
<b>Cy3</b>	X	X	-
<b>ATTO 532</b>	X	-	-
<b>5/6-Rhodamine</b>	X	-	X
<b>ATTO 488</b>	X	-	X
<b>5/6-Fluorescein</b>	X	-	X
<b>AF488</b>	X	-	X
<b>ATTO 425</b>	X	-	X
<b>DEAC</b>	X	-	X
<b>AMCA</b>	-	-	-
<b>AF405</b>	-	-	-