



Click Chemistry describes pairs of functional groups that rapidly react (“click”) with each other under mild, aqueous conditions. **Molecules** that have been **functionalized with a CLICK moiety** (e.g. a terminal Alkyne) **can be detected by labeling with a compatible CLICKable fluorescent dye** (e.g. a fluorescent Azide) that allows the **subsequent visualization by fluorescence spectroscopy**.

A number of CLICKable fluorescent dyes have been thoroughly selected to cover the whole UV-VIS spectrum:

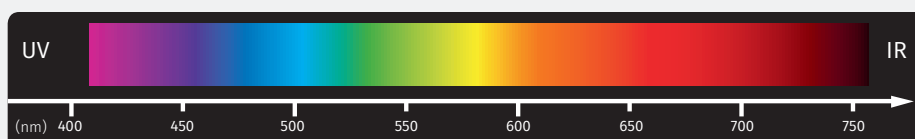


Figure 1: Visible spectrum of light.

Go to www.jenabioscience.com/clickabledyes

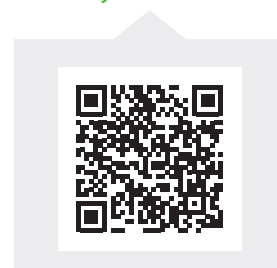


Table 1: Overview of available CLICKable fluorescent dyes. | ✓ = Product link. Please click for more information.

Emission Color	Labeling dye	Azide	Picolyl-Azide	Alkyne	DBCO* =ADIBO/DIBAC	Tetrazine	6-Methyl-Tetrazine
blue	7-Hydroxycoumarin	✓					
	ATTO425				✓		
green	5-Fluorescein	✓		✓		✓	✓
	6-Fluorescein	✓		✓			
	5/6-Fluorescein		✓		✓		
	Dansyl	✓					
	5/6-Carboxyrh. 110	✓		✓	✓		
	ATTO 488				✓	✓	✓
	AF488 (also known as Alexa Fluor®488)	✓	✓	✓	✓		
	BDP-FL (also known as BODIPY®FL)	✓		✓	✓	✓	✓
yellow green	ATTO 532					✓	✓
yellow	Sulfo-Cy3	✓	✓	✓	✓		✓
	Cy3	✓				✓	
	5-TAMRA	✓		✓		✓	✓
	5/6-TAMRA	✓	✓	✓	✓		
	5/6-Sulforhodamine B	✓		✓	✓		
orange	5/6-Texas Red	✓		✓	✓		
	AF594 (also known as Alexa Fluor®594)	✓	✓	✓	✓		
red	Sulfo-Cy5	✓	✓	✓	✓		✓
	Cy5	✓				✓	
	ATTO 647N					✓	✓
	Cy5.5	✓	✓	✓	✓		
Near-IR	Cy7	✓	✓	✓	✓		

*DBCO = Dibenzocyclooctyne, ADIBO = Azadibenzocyclooctyne, DIBAC = Dibenzozacyclooctyne

Check out our complete Click Chemistry product portfolio and find more information at www.click-chemistry.net

