

Product Information

Cyanine Dye, Alkyne

Unit Size: 1 mg

Technical Summary

Catalog no.	Dye	Ex/Em(nm)	MW
92100	Cyanine 555	555/565	769
92101	Cyanine 647	650/665	795

Storage and Handling

Store cyanine dye, alkyne at -20°C, protected from moisture and light. Product is stable for at least 12 months from date of receipt if stored as recommended.

Stock solutions may be prepared in water, DMF, or DMSO. For making stock solutions, we recommend dissolving the dye in water at 10 mM. Stock solutions prepared in water can be stored at \leq -20°C for at least 12 months.

Product Description

Red fluorescent Cyanine 555 alkyne and far-red fluorescent Cyanine 647 alkyne react with azide to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition. Both Cyanine 555 alkyne and Cyanine 647 alkyne have been used to analyze O-GlcNAcylated proteins in "2D glyco-DIGE" platform.

Figure 1. Cyanine 555 Alkyne

Figure 2. Cyanine 647 Alkyne

Related Products

Catalog number	Product	
90082	DMSO, Anhydrous	
90016 90049	Cyanine Dye, Succinimidyl Ester	
92086 96006	CF® Dye Alkyne	
92103 96095	CF® Dye SE/TFP Ester	
92230 92450	Mix-n-Stain™ CF® Dye Antibody Labeling Kits	
22014	Ultrafiltration Vial, 10K MWCO (5 per pack)	
22018	Ultrafiltration Vial, 3K MWCO (5 per pack)	

Please visit www.biotium.com to view our full selection of CF® reactive dyes and labeling kits, CF® Dye labeled antibodies and other conjugates, and more.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.