Cyanine Dye Alkyne

Cyanine 555 and Cyanine 647 alkynes can be used for the analysis of O-GlcNAcylated proteins by 2D glyco-DIGE.



Product attributes

Call us: 800-304-5357

Chemical reactivity (reacts with)	Azides/Picolyl azides
Functional group	Alkyne
Storage Conditions	Store at -10 to -35 °C, Protect from light

Email: btinfo@biotium.com

Product Description

Cyanine 555 and Cyanine 647 alkynes can be used to analyze O-GlcNAcylated proteins by 2D glyco-DIGE (differential gel electrophoresis). They react with azide groups to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition.

- Size- and charge-matched for analysis of O-GlcNAcylated proteins via 2D glyco-DIGE.
- React with azides to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition.

Cyanine 555:

- $\lambda_{Ex}/\lambda_{Em}$ (water) = 555/565 nm
 - Dark red solid soluble in water, DMF and DMSO
 - C40H56N4O7S2
 - MW: 769

Cyanine 647:

- λEx/λEm (water) = 650/665 nm
 - Dark blue solid soluble in water, DMF and DMSO
 - C42H58N4O7S2
 - MW: 795

Also see CF® Dye Alkynes, available with a wide selection of our bright and photostable CF® dyes.

This datasheet was generated on November 4, 2025 at 12:56:40 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cyanine-dye-alkyne/