

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

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

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
- $C_{40}H_{56}N_4O_7S_2$
- MW: 769

Cyanine 647:

- $\lambda_{Ex}/\lambda_{Em}$ (water) = 650/665 nm
- Dark blue solid soluble in water, DMF and DMSO
- $C_{42}H_{58}N_4O_7S_2$
- MW: 795

Also see [CF® Dye Alkynes](#), available with a wide selection of our bright and photostable CF® dyes.