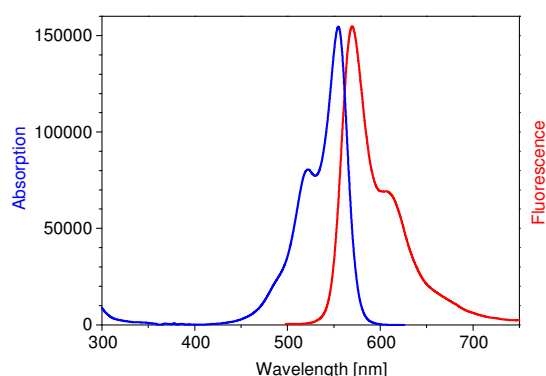


DBCO-SETA 555

Cat. No.	Amount
CLK-A126-05	0,5 mg



Absorption and emission spectrum of DBCO-SETA 555 in phosphate buffer (pH 7.4)

Molecular Weight: 1147.32 g/mol (protonated form)

Spectroscopic properties:

λ_{abs} 555 nm; λ_{em} 570 nm; ϵ 155.000 cm⁻¹ M⁻¹;
 quantum yield 7 % (in phosphate buffer pH 7.4)

Storage conditions: store at -20 °C and dark

Shelf life: 12 months

Purity: > 90 % (HPLC)

Appearance: dark red solid

Solubility: DMSO, DMF, methanol, water

Insoluble: acetone, chloroform, toluene

For research use only!

Product Features and Benefits:

DBCO-SETA 555 is a highly hydrophilic click chemistry reagent containing one dibenzocyclooctyne (DBCO) group for strain-mediated cycloaddition reactions (copper-free click reactions) with azide-modified reagents and biomolecules. It features the same excitation and emission wavelengths as Cy3, Alexa Fluor™ 555 or Seta 555 and can therefore be used with these filter sets. It combines high photostability and brightness.

Advantages

- Perfectly suited for excitation with the 532-nm laser
- Good aqueous solubility: this label does not alter the solubility of the protein conjugate
- Low molecular weight: DBCO-SETA 555 does not add substantial mass to the conjugates
- Photostability: Higher photostability as compared to Alexa Fluor™ 555 or Cy3
- Ideal for non-radioactive labeling of azide modified molecules (drugs, proteins, oligos and DNA)