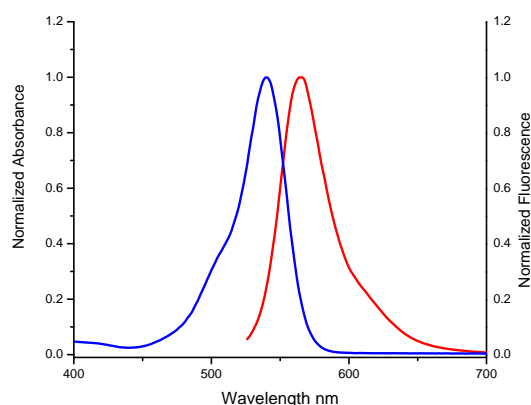
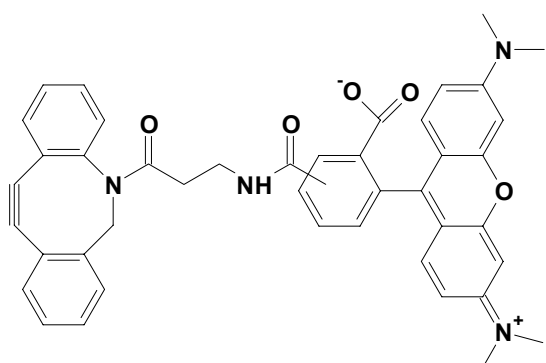


## Dibenzylcyclooctyne-Fluor 545

Cat. No.	Amount
CLK-A110-2	2 mg
CLK-A110-5	5 mg



Absorption and emission spectrum Fluor 545

**Molecular formula:** C<sub>43</sub>H<sub>36</sub>N<sub>4</sub>O<sub>5</sub>

**Molecular weight:** 688.78 g/mol

**Spectroscopic properties:**

$\lambda_{\text{abs}}$  546 nm;  $\lambda_{\text{em}}$  565 nm,  $\epsilon$  92000 cm<sup>-1</sup> M<sup>-1</sup> (in MeOH)

**Storage conditions:** store at -20°C

**Purity:** >90% (H NMR)

**Appearance:** dark red solid

**Shelf life:** 12 months

**Solubility:** DMSO, DMF, DCM, Chloroform, MeOH

**Product Features and Benefits:**

With excitation maximum at 546 nm, the dibenzylcyclooctyne – conjugate is an excellent match to the intense 546 nm spectral line of the mercury-arc lamps used in most fluorescence microscopes. This probe is also efficiently excited by the 543 nm spectral line of the Ar – Kr mixed gas laser commonly used in many confocal laser-scanning microscopes.

The fluorescence of the dibenzylcyclooctyne – tetramethylrhodamine conjugate as well as of many other rhodamine based probes is pH insensitive between 4 and 9, and exhibits excellent photostability.

**For research use only!**