

5-IAF [5-lodoacetamidofluorescein] *CAS 63368-54-7*

Catalog number: 222 Unit size: 25 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Brown solid

Molecular Weight 515.25

Soluble In DMSO

Chemical Structure

Spectral Properties

Excitation Wavelength 498 nm

Emission Wavelength 517 nm

Applications

5-IAF is one of the most popular green fluorescent dyes for thiol modifications of proteins along with fluorescein-5-maleimide. The resulted bioconjugates are quite stable.