

Product Information Sheet

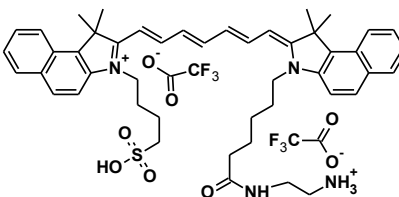
Ordering Information

Product Number: 188
Product Name: ICG amine
Unit Size: 1 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 24 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 1001.08

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 780 nm
Emission Wavelength: 800 nm

Application Notes

Indocyanine green (ICG) is a cyanine dye used in medical diagnostics. It is used for determining cardiac output, hepatic function, and liver blood flow, and for ophthalmic angiography. It has a peak spectral absorption close to 800 nm. These infrared frequencies penetrate retinal layers, allowing ICG angiography to image deeper patterns of circulation than fluorescein angiography. This ICG amine can be used to modify the molecules that contain a carbonyl group (such as an aldehyde or a carboxy group).