

ICG acid

Catalog number: 189 Unit size: 5 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Green solid
Molecular Weight	844.98
Soluble In	DMSO
Chemical Structure	
Spectral Properties	

Spectral Properties

Applications

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 acid can be used to modify the molecules that contain an amino group.