

# GTIF488 [iFluor® 488 gentamicin conjugate]

Catalog Number: 24302, 24303

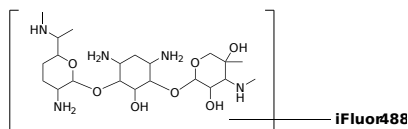
Unit Size: 100 ug, 1 mg

## Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

Appearance	Solid red
Molecular Weight	~1100
Soluble In	DMSO
Chemical Structure	



## Spectral Properties

Excitation Wavelength	491 nm
Emission Wavelength	516 nm

## Applications

Gentamicin is widely used for treating tuberculosis and Gram-negative infections and is particularly useful in neonatal intensive care units. Fluorophore-tagged drugs, such as gentamicin- iFluor® 488 (GTIF488) conjugates, might be used to identify drug trafficking routes across the BLB and into hair cells, and specifically in the cochlea - the sensory organ responsible for hearing. Identifying the mechanisms involved in the intracochlear trafficking of ototoxic drugs to hair cells is of fundamental as well as clinical importance. GTIF488 is well excited with the common Argon laser at 488 nm. It provides researchers a valuable tool to study the cellular uptake, kinetics and distribution of gentamicin through fluorescence techniques, such as fluorescence microscopy, fluorescence correlation spectroscopy, single molecule spectroscopy. The iFluor® 488 is well compatible with the common FITC filter set.