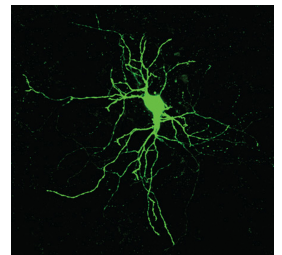
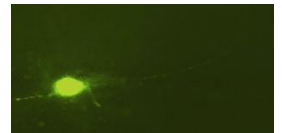
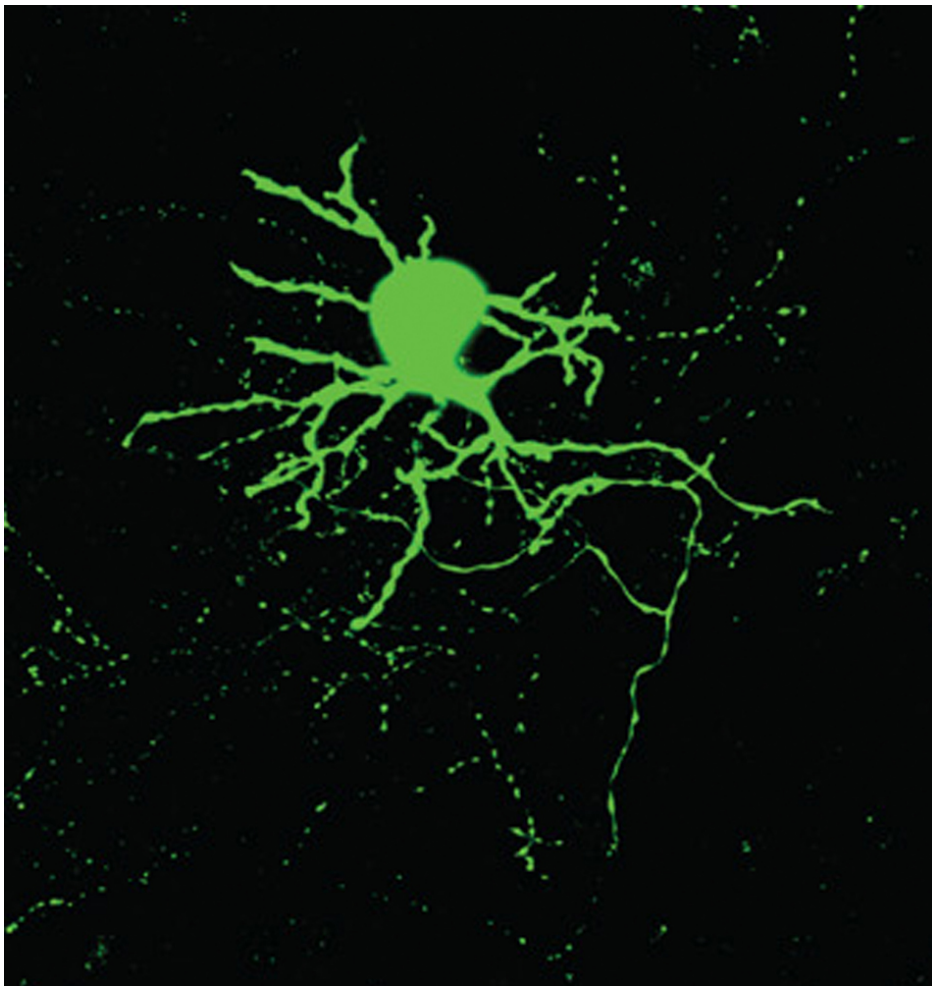


# NEUROBIOTIN® 488 Tracer

## Product Images



## Short Description

---

NEUROBIOTIN 488 Tracer is a tri-functional molecule designed for neuronal tracing and cell filling.

### Features:

- Bright green fluorophore, similar in fluorescence to fluorescein, Cy2 or Alexa Fluor® 488
- Biotin label with a biotinidase-resistant linkage
- Fixable primary amine
- Used for visualizing neural architecture and for the identification of gap junction coupling
- Can be used in many types of preparations including in vivo, whole mounts, slice preparations, or cultured cells
- Can be delivered by many routes such as intracellular electrodes, microinjection, cut-loading, or scrape-loading

### Advantages of NEUROBIOTIN Tracer over biocytin and other neuronal labels:

- Better solubility
- More efficiently iontophoresed
- Remains in cell longer
- Non-toxic
- Can be fixed with formalin or glutaraldehyde

# Additional Information

Unit Size	2 mg
Maximum Emission	517 nm
Recommended Storage	2-8 °C (desiccated). Once in solution, store frozen. This product does not contain an antimicrobial agent.
Recommended Usage	NEUROBIOTIN 488 is soluble under the following conditions.Solvents and concentrations will require optimization depending on the application. Up to 10% in 3.0 M Tris, pH 8.8 Up to 2% in 0.25 M Tris, pH 8.8 or 0.25 M Tris, 1.0 M NaCl, pH 8.8 Up to 2% in 0.1 M bicarbonate buffer, pH 8.5 or 0.25 M bicarbonate, 1.0 M NaCl, pH 8.5 Up to 0.5% in 0.13 M cesium gluconate, 10 mM HEPES, pH 7.2 Up to 10% in DMSO
Molecular Weight	747
Maximum Excitation	493 nm
Neuronal Tracer - Direction of Transport	Anterograde/Retrograde
Detection Method	Avidin(Streptavidin)/Biotin Method, Chromogenic, Fluorescence

