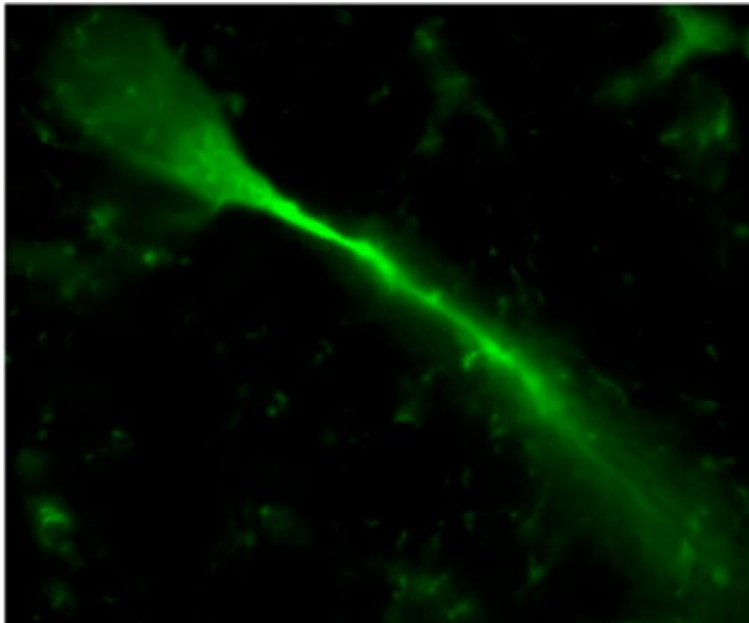




# NEUROBIOTIN®-PLUS Tracer

## Product Images



## Short Description

NEUROBIOTIN-Plus Tracer is a biotin label that is impervious to cleavage and breakdown by biotinidase. It is used for visualizing neural architecture and for the identification of gap junction coupling

### Features:

- Fixable primary amine
- 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
- Biotin label can be detected using avidin or streptavidin systems with either chromogenic or fluorescence visualization methods

### 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	5 mg
Recommended Storage	2-8 °C (desiccated). Once in solution, store frozen. This product does not contain an antimicrobial agent.
Recommended Usage	Fixable with formaldehyde or glutaraldehyde
Solution	10% solutions can be made in water, buffers, 1 M KCl, 1 M NaCl, or dimethylsulfoxide.
Molecular Weight	358
Neuronal Tracer - Direction of Transport	Anterograde/Retrograde
Detection Method	Avidin(Streptavidin)/Biotin Method, Chromogenic, Fluorescence

