

PRODUCT INFORMATION



iFluor™ 800-Streptavidin Conjugate

Item No. 44616

Ex./Em. Max: 801/820 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Light sensitive

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

iFluor™ 800-streptavidin conjugate is a fluorescent conjugate composed of iFluor™ 800 conjugated to streptavidin. It displays excitation/emission maxima of 801/820 nm, respectively. iFluor™ 800-streptavidin conjugate can be used as a second-step reagent for indirect immunofluorescent staining when used in conjunction with a biotinylated primary antibody.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/24/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM