

PRODUCT INFORMATION



iFluor® 488 Azide

Item No. 44066

FW: 716.6
Ex./Em. Max.: 491/516 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.

Laboratory Procedures

iFluor® 488 azide is supplied as a solid. A stock solution may be made by dissolving the iFluor® 488 azide in the solvent of choice, which should be purged with an inert gas. iFluor® 488 azide is soluble in DMSO.

Description

iFluor® 488 azide is a fluorophore functionalized with azide, which reacts with alkynes, such as dibenzocyclooctyne (DBCO), to form a stable conjugate. It displays excitation/emission maxima of 491/516 nm, respectively. iFluor® 488 azide can be used for labeling alkyne-modified biological molecules for fluorescence imaging and other fluorescence-based biological applications. It has been used to detect 5-ethynyl-2'-deoxyuridine (EdU; Item No. 20518) incorporation into newly synthesized DNA for the quantification of cell proliferation.¹

Reference

1. Chang, C.-J. and Taniguchi, A. Establishment of a nanopatterned renal disease model by mimicking the physical and chemical cues of a diseased mesangial cell microenvironment. *ACS Appl. Bio Mater.* **4**(2), 1573-1583 (2021).

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, 11/25/2025

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