

PRODUCT INFORMATION



iFluor® 647 TCO

Item No. 44074

Synonym: iFluor® 647 *trans*-cyclooctene

FW: 1,181.3

Purity: ≥90%

Ex./Em. Max: 656/670 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® 647 TCO is supplied as a solid. A stock solution may be made by dissolving the iFluor® 647 TCO in the solvent of choice, which should be purged with an inert gas. iFluor® 647 TCO is soluble in DMSO.

Description

iFluor® 647 TCO is a fluorophore functionalized with *trans*-cyclooctene (TCO), which reacts with tetrazine rings to form a stable conjugate. It displays excitation/emission maxima of 656/670 nm, respectively. iFluor® 647 TCO can be used for labeling tetrazine-modified biological molecules for fluorescence imaging and other fluorescence-based biological applications.

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, 04/23/2026

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