

PRODUCT INFORMATION



LDL-DyLight® 488 Assay Reagent

Item No. 601471

Synonyms:	Low-density Lipoprotein-DyLight® 488
Contents:	This vial contains 120 µl of human LDL conjugated to DyLight® 488 in PBS. DyLight® 488 is a product of Thermo Fisher Scientific Inc.
Ex./Em. Max:	488/518 nm
Storage:	4°C
Stability:	≥18 months

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

Description

LDL-DyLight® 488 Assay Reagent is composed of LDL conjugated to DyLight® 488. It displays excitation/emission maxima of 488/518 nm, respectively, and can be used to visualize the uptake of LDL in live cells. LDL-DyLight® 488 Assay Reagent is also available as part of Cayman's LDL Uptake Flow Cytometry Assay Kit (Item No. 601470).

Instructions for Use

Mix the tube of LDL-DyLight® 488 Assay Reagent gently (*do not vortex!*) and dilute 1:10-1:500 in cell culture medium.

NOTE: If particulates are observed, the diluted reagent can be filtered using a 0.45 µm syringe-top filter.

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, 12/09/2025

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