

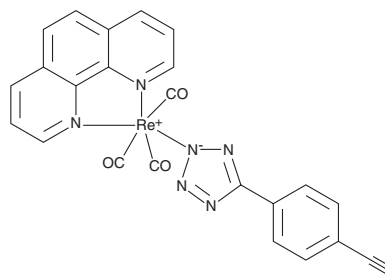
PRODUCT INFORMATION



ReZolve-Alkyne™

Item No. 25909

CAS Registry No.: 2070860-85-2
Formal Name: (OC-6-33)-tricarbonyl[5-(4-ethynylphenyl)-2H-tetrazolato-κN²](1,10-phenanthroline-κN¹,κN¹⁰)-rhenium
MF: C₂₄H₁₃N₆O₃Re
FW: 619.6
Ex./Em. Max: 405/570 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Description

ReZolve-Alkyne™ is a fluorogenic probe that can be used to tag azide-functionalized molecules in cells and tissues. The terminal alkyne on ReZolve-Alkyne™ can be coupled to molecules of interest using click chemistry in the presence of a copper catalyst and visualized using epifluorescent, confocal, and two-photon microscopy. It displays excitation/emission maxima of 405/570 nm, respectively.

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, 07/26/2018

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