

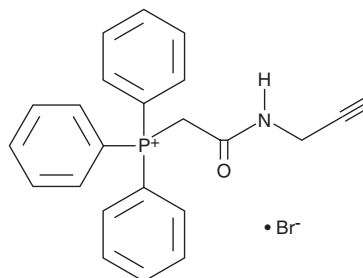
PRODUCT INFORMATION



WYneN

Item No. 42010

CAS Registry No.: 3052492-38-0
Formal Name: (2-oxo-2-(prop-2-yn-1-ylamino)ethyl) triphenylphosphonium, monobromide
Synonym: Wittig-alkyne N
MF: C₂₃H₂₁NOP • Br
FW: 438.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

WYneN is supplied as a solid. A stock solution may be made by dissolving the WYneN in the solvent of choice, which should be purged with an inert gas. WYneN is sparingly soluble (1-10 mg/ml) in the organic solvent DMSO.

Description

WYneN is a chemoselective probe for the detection of sulfenylated proteins in live cells.¹ It is selective for sulfenylated proteins over 5,5'-dithiobis-(2-nitrobenzoic acid) (DTNB) and 4,4'-dipyridyl disulfide (4-DPS). WYneN binds to sulfenylated proteins in live cells and can then be conjugated to azide-containing fluorescent probes for detection.

Reference

1. Shi, Y., Fu, L., Yang, J., *et al.* Wittig reagents for chemoselective sulfenic acid ligation enables global site stoichiometry analysis and redox-controlled mitochondrial targeting. *Nat. Chem.* **13(11)**, 1140-1150 (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, 10/25/2024

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