

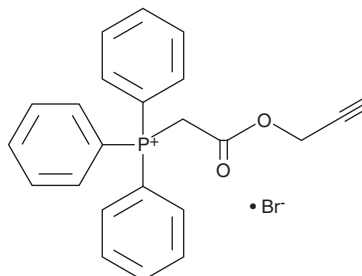
PRODUCT INFORMATION



WYneO

Item No. 42011

CAS Registry No.: 1648555-11-6
Formal Name: [2-oxo-2-(2-propyn-1-yloxy)ethyl] triphenyl-phosphonium, monobromide
Synonym: Wittig-alkyne O
MF: $C_{23}H_{20}O_2P \cdot Br$
FW: 439.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of WYneO can be prepared by directly dissolving the solid in aqueous buffers. WYneO is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

WYneO is a probe for the detection of sulfenylated proteins.^{1,2}

References

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).
2. He, J.-X., Fei, Z.-C., Fu, L., *et al.* A modification-centric assessment tool for the performance of chemoproteomic probes. *Nat. Chem. Biol.* **18**(8), 904-912 (2022).

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, 10/18/2024

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