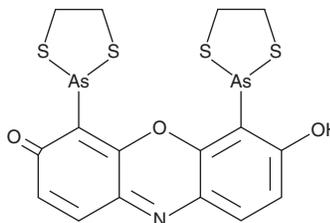


PRODUCT INFORMATION



ReAsH-EDT₂ Item No. 19767

CAS Registry No.: 438226-89-2
Formal Name: 4,6-bis(1,3,2-dithiarsolan-2-yl)-7-hydroxy-3H-phenoxazin-3-one
Synonym: Lumio Red
MF: C₁₆H₁₃As₂NO₃S₄
FW: 545.4
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

ReAsH-EDT₂ is supplied as a solid. A stock solution may be made by dissolving the ReAsH-EDT₂ in the solvent of choice, which should be purged with an inert gas. ReAsH-EDT₂ is slightly soluble in chloroform.

Description

ReAsH-EDT₂ is a fluorescent, membrane-permeable biarsenical compound that binds covalently to tetracysteine sequences, which are engineered into target proteins.¹ It binds proteins that have the tetracysteine tag almost immediately after translation.¹ ReAsH-EDT₂ is commonly used to study protein trafficking, folding, and interactions in living cells or cell lysates.¹⁻³ This red-emitting fluorophore is excited at 593 nm, with emission at 608 nm.³

References

1. Rudner, L., Nydegger, S., Coren, L. V., *et al.* Dynamic fluorescent imaging of human immunodeficiency virus type 1 gag in live cells by biarsenical labeling. *J. Virol.* **79**(7), 4055-4065 (2005).
2. Luedtke, N. W., Dexter, R. J., Fried, D. B., *et al.* Surveying polypeptide and protein domain conformation and association with FIAsH and ReAsH. *Nat. Chem. Biol.* **3**(12), 779-784 (2007).
3. Perdios, L., Bunney, T. D., Warren, S. C., *et al.* Time-resolved FRET reports FGFR₁ dimerization and formation of a complex with its effector PLCγ1. *Adv. Biol. Regul.* **60**, 6-13 (2016).

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, 09/29/2022

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