

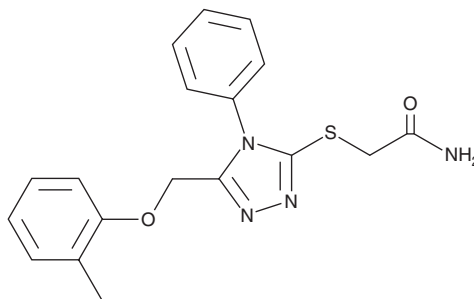
PRODUCT INFORMATION



A2ti-2

Item No. 43719

CAS Registry No.: 482646-13-9
Formal Name: 2-[[5-[(2-methylphenoxy)methyl]-4-phenyl-4H-1,2,4-triazol-3-yl]thio]-acetamide
MF: C₁₈H₁₈N₄O₂S
FW: 354.4
Purity: ≥98%
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

A2ti-2 is supplied as a solid. A stock solution may be made by dissolving the A2ti-2 in the solvent of choice, which should be purged with an inert gas. A2ti-2 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

A2ti-2 is an inhibitor of the protein-protein interaction between annexin A2 and S100A10 (IC₅₀ = 230 μM in a cell-free assay).¹ It inhibits human papillomavirus (HPV) entry in HeLa cervical cancer cells when used at concentrations of 50, 75, or 100 μM.²

References

1. Reddy, T.R.K., Li, C., Fischer, P.M., *et al.* Three-dimensional pharmacophore design and biochemical screening identifies substituted 1,2,4-triazoles as inhibitors of the annexin A2-S100A10 protein interaction. *ChemMedChem* **7**(8), 1435-1446 (2012).
2. Woodham, A.W., Taylor, J.R., Jimenez, A.I., *et al.* Small molecule inhibitors of the annexin A2 heterotetramer prevent human papillomavirus type 16 infection. *J. Antimicrob. Chemother.* **70**(6), 1686-1690 (2015).

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, 10/08/2025

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