

PRODUCT INFORMATION



Maleimide-PEG₂-AZD

Item No. 20129

CAS Registry No.: 1037589-69-7
Formal Name: 2,5-dihydro-2,5-dioxo-N-[2-[2-[3-oxo-3-[[4-[3-oxo-3-(2-oxo-1-azetidiny)]propyl]phenyl]amino]propoxy]ethoxy]ethyl]-1H-pyrrole-1-propanamide

Synonym: Mal-PEG₂-AZD

MF: C₂₆H₃₂N₄O₈

FW: 528.6

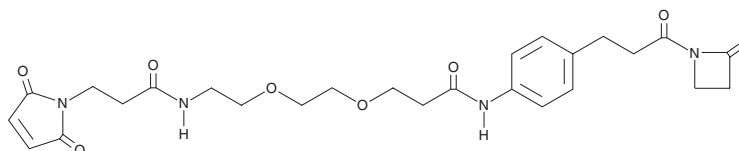
Purity: ≥98%

UV/Vis.: λ_{max}: 248 nm

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

Maleimide-PEG₂-AZD is supplied as a solid. A stock solution may be made by dissolving the maleimide-PEG₂-AZD in the solvent of choice, which should be purged with an inert gas. Maleimide-PEG₂-AZD is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of maleimide-PEG₂-AZD in these solvents is approximately 30 mg/ml.

Maleimide-PEG₂-AZD is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, maleimide-PEG₂-AZD should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Maleimide-PEG₂-AZD has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Maleimide-PEG₂-AZD is a heterobifunctional bioconjugation linker.^{1,2} It has been used in the development of protein- and peptidomimetic-antibody conjugates, including the FGF21 protein-antibody conjugate PF-05231023 and κ-opioid receptor (KOR) agonists, respectively.

References

1. Roberts, L.R., Brady, K.D., Brown, A., *et al.* Kappa agonist CovX-Bodies. *Bioorg. Med. Chem. Lett.* **22**(12), 4173-4178 (2012).
2. Dirksen, A., Davis, K.A., Collins, J.T., *et al.* Process development of a FGF21 protein-antibody conjugate. *Peptide Sci.* **110**(1), e23042 (2017).

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, 04/11/2025

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