

PRODUCT INFORMATION



UK 383367

Item No. 15029

CAS Registry No.: 348622-88-8

Formal Name: (βR)-3-(aminocarbonyl)-β-(3-cyclohexylpropyl)-N-hydroxy-1,2,4-oxadiazole-5-propanamide

MF: C₁₅H₂₄N₄O₄

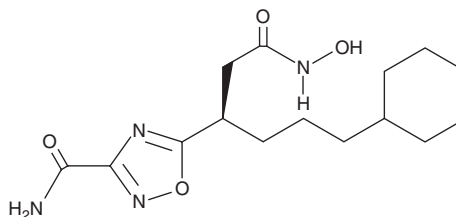
FW: 324.4

Purity: ≥98%

Supplied as: A crystalline 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

UK 383367 is supplied as a crystalline solid. A stock solution may be made by dissolving the UK 383367 in the solvent of choice, which should be purged with an inert gas. UK 383367 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UK 383367 in ethanol and DMF is approximately 16 mg/ml and approximately 25 mg/ml in DMSO.

UK 383367 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UK 383367 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. UK 383367 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UK 383367 is a potent inhibitor of bone morphogenetic protein 1 (BMP1; IC₅₀ = 44 nM).^{1,2} It is more than 200-fold selective for BMP1 over a panel of matrix metalloproteinases. UK 383367 is effective in a cell-based model of collagen deposition and is very effective at penetrating human skin *in vitro*.²

References

1. Fish, P. V., Allan, G. A., Bailey, S., *et al.* Potent and selective nonpeptidic inhibitors of procollagen C-proteinase. *J. Med. Chem.* **509(15)**, 3442-3456 (2007).
2. Bailey, S., Fish, P. V., Billotte, S., *et al.* Succinyl hydroxamates as potent and selective non-peptidic inhibitors of procollagen C-proteinase: Design, synthesis, and evaluation as topically applied, dermal anti-scarring agents. *Bioorg. Med. Chem. Lett.* **18(24)**, 6562-6567 (2008).

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, 12/06/2022

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