

PRODUCT INFORMATION

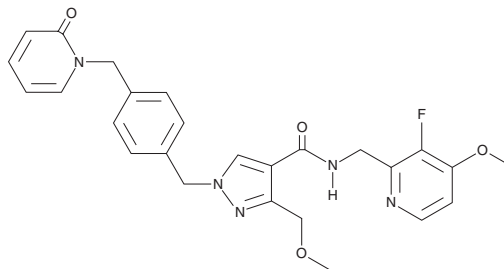


Sebetralstat

Item No. 45043

CAS Registry No.: 1933514-13-6
Formal Name: N-[(3-fluoro-4-methoxy-2-pyridinyl)methyl]-3-(methoxymethyl)-1-[[4-[(2-oxo-1(2H)-pyridinyl)methyl]phenyl]methyl]-1H-pyrazole-4-carboxamide

Synonym: KVD900
MF: C₂₆H₂₆FN₅O₄
FW: 491.5
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

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

Description

Sebetralstat is an inhibitor of plasma kallikrein (IC₅₀ = 6 nM).¹ It is selective for plasma kallikrein over a panel of 14 additional proteases (IC₅₀s = >10 μM for all). Sebetralstat (50-1,000 nM) inhibits dextran sulfate-induced cleavage of kininogen-1 in isolated human plasma. Formulations containing sebetralstat have been used in the treatment of acute hereditary angioedema (HAE).

Reference

1. Duckworth, E.J., Murugesan, N., Li, L., *et al.* Pharmacological suppression of the kallikrein kinin system with KVD900: An orally available plasma kallikrein inhibitor for the on-demand treatment of hereditary angioedema. *Clin. Exp. Allergy* **52(9)**, 1059-1070 (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, 02/06/2026

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