

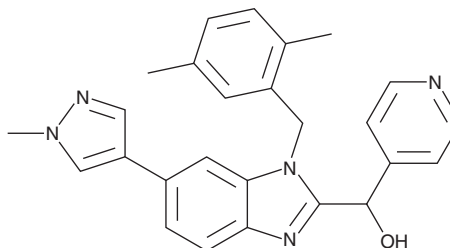
PRODUCT INFORMATION



UCB 9260

Item No. 37336

CAS Registry No.: 1515888-53-5
Formal Name: 1-[(2,5-dimethylphenyl)methyl]-6-(1-methyl-1H-pyrazol-4-yl)- α -4-pyridinyl-1H-benzimidazole-2-methanol
MF: C₂₆H₂₅N₅O
FW: 423.5
Purity: $\geq 95\%$
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

UCB 9260 is supplied as a solid. A stock solution may be made by dissolving the UCB 9260 in the solvent of choice, which should be purged with an inert gas. UCB 9260 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of UCB 9260 in these solvents is approximately 20 mg/ml. UCB 9260 is slightly soluble in ethanol.

Description

UCB 9260 is a TNF inhibitor.¹ It binds to trimeric TNF ($K_d = 13.8$ nM) and selectively increases the thermal stability of the TNF trimer over 10 other TNF superfamily members at 100 μ M. UCB 9260 inhibits TNF-induced NF- κ B activity in a reporter assay using HEK293 cells and TNF-induced cytotoxicity in L-929 fibroblasts (IC_{50} s = 202 and 116 nM, respectively). *In vivo*, UCB 9260 (30-300 mg/kg, p.o.) reduces TNF-induced peritoneal neutrophil recruitment in mice. Intraperitoneal administration of UCB 9260 (150 mg/kg) decreases disease severity in a mouse model of collagen antibody-induced arthritis.

Reference

1. O'Connell, J., Porter, J., Kroepfli, B., *et al.* Small molecules that inhibit TNF signalling by stabilising an asymmetric form of the trimer. *Nat. Commun.* **10**(1), 5795 (2019).

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, 8/22/2023

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