

PRODUCT INFORMATION

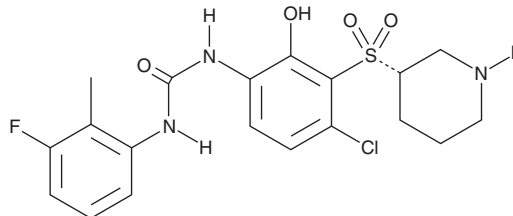


Danirixin

Item No. 20980

CAS Registry No.: 954126-98-8
Formal Name: N-[4-chloro-2-hydroxy-3-[(3S)-3-piperidinylsulfonyl]phenyl]-N'-(3-fluoro-2-methylphenyl)-urea

Synonym: GSK1325756
MF: C₁₉H₂₁ClFN₃O₄S
FW: 441.9
Purity: ≥98%
UV/Vis.: λ_{max}: 266, 323 nm
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.

Description

Danirixin is an antagonist of chemokine (C-X-C motif) receptor 2 (CXCR2; IC₅₀ = 12.5 μM).¹ It is selective for CXCR2 over CXCR1 (IC₅₀ = 977 nM). Danirixin inhibits IL-8-induced calcium mobilization in CHO-K1 cells expressing human CXCR2 (pA₂ = 7.52). It inhibits neutrophil surface expression of CD11b in isolated rat whole blood (IC₅₀ = 890 nM). Danirixin reduces bronchoalveolar lavage fluid (BALF) neutrophil infiltration in rat models of LPS- or ozone-induced pulmonary neutrophilia (ED₅₀s = 1.4 and 16 mg/kg, respectively).

Reference

1. Busch-Petersen, J., Carpenter, D.C., Burman, M., *et al.* Danirixin: A reversible and selective antagonist of the CXC chemokine receptor 2. *J. Pharmacol. Exp. Ther.* **362**(2), 338-346 (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, 10/19/2022

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