

PRODUCT INFORMATION

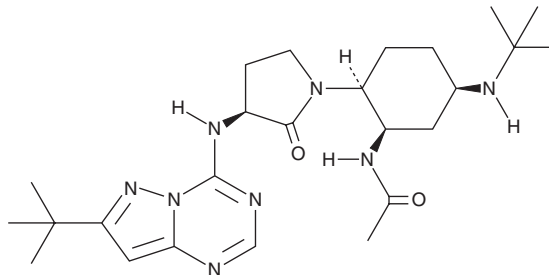


BMS 813160

Item No. 28421

CAS Registry No.: 1286279-29-5
Formal Name: N-[(1R,2S,5R)-5-[(1,1-dimethylethyl)amino]-2-[(3S)-3-[[7-(1,1-dimethylethyl)pyrazolo[1,5-a]-1,3,5-triazin-4-yl]amino]-2-oxo-1-pyrrolidinyl]cyclohexyl]-acetamide

MF: C₂₅H₄₀N₈O₂
FW: 484.6
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 272 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.

Laboratory Procedures

BMS 813160 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 813160 in the solvent of choice, which should be purged with an inert gas. BMS 813160 is soluble in the organic solvent DMSO.

Description

BMS 813160 is a dual antagonist of chemokine receptor 2 (CCR2) and CCR5 (IC₅₀s = 6.2 and 3.6 nM, respectively, in radioligand binding assays).^{1,2} It inhibits chemotaxis induced by the CCR2 ligand MCP-1 in THP-1 cells the CCR5 ligand MIP-1β in isolated peripheral T cells with IC₅₀ values of 0.8 and 1.1 nM, respectively.¹ BMS 813160 inhibits peritoneal monocyte and macrophage infiltration in an hCCR-2 KI mouse model of thioglycolate-induced peritonitis, with 54% inhibition when administered at a dose of 50 mg/kg twice per day.

References

1. Carter, P.H., Cherney, R.J., Rosso, V.W., *et al.* N-((1r,2s,5r)-5-(tert-butylamino)-2-((s)-3-(7-tert-butylpyrazolo[1,5-a][1,3,5]triazin-4-ylamino)-2-oxopyrrolidin-1-yl)cyclohexyl)acetamide, a dual modulator of chemokine receptor activity, crystalline forms and processes. **PCT /US20 I 0/0 52288** (2011).
2. Norman, P. A dual CCR2/CCR5 chemokine antagonist, BMS-813160? Evaluation of WO2011046916. *Expert. Opin. Ther. Pat.* **21(12)**, 1919-1924 (2011).

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, 12/05/2022

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