

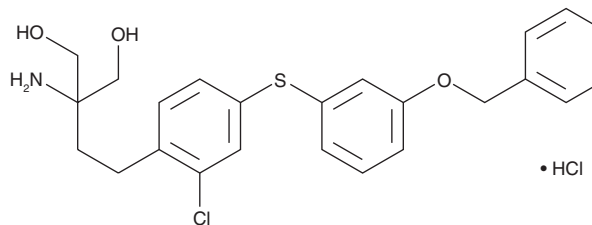
PRODUCT INFORMATION



KRP 203

Item No. 10010426

CAS Registry No.: 509088-69-1
Formal Name: 2-amino-2-[2-[2-chloro-4-[[3-(phenylmethoxy)phenyl]thio]phenyl]ethyl]-1,3-propanediol, monohydrochloride
Synonym: Mocravimod (hydrochloride)
MF: C₂₄H₂₆ClNO₃S • HCl
FW: 480.4
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

KRP 203 is supplied as a crystalline solid. A stock solution may be made by dissolving the KRP 203 in an organic solvent purged with an inert gas. KRP 203 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of KRP 203 in these solvents is approximately 15 and 30 mg/ml, respectively.

KRP 203 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KRP 203 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. KRP 203 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KRP 203 is a sphingosine-1-phosphate receptor 1 (S1P₁) agonist.¹ It selectively induces calcium mobilization in CHO-K1 cells expressing mouse S1P₁ over mouse S1P₃ and S1P₄ (EC₅₀s = 0.84, >1,000, and 9.61 nM, respectively). *In vivo*, KRP 203 (0.3 mg/kg) prevents the development of wasting syndrome, decreases diarrhea and colonic inflammatory cell infiltration, and increases colon length in an *IL10*^{-/-} mouse model of colitis. KRP 203 inhibits inflammatory cell infiltration and prolongs allograft survival in a rat model of minor histocompatibility complex-disparate heart transplantation.²

References

1. Song, J., Matsuda, C., Kai, Y., *et al.* A novel sphingosine 1-phosphate receptor agonist, 2-amino-2-propanediol hydrochloride (KRP-203), regulates chronic colitis in interleukin-10 gene-deficient mice. *J. Pharmacol. Exp. Ther.* **324**(1), 276-283 (2007).
2. Shimizu, H., Takahashi, M., Kaneko, T., *et al.* KRP-203, a novel synthetic immunosuppressant, prolongs graft survival and attenuates chronic rejection in rat skin and heart allografts. *Circulation* **111**, 222-229 (2005).

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, 08/11/2023

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