

PRODUCT INFORMATION



Rimiducid

Item No. 35303

CAS Registry No.: 195514-63-7

Formal Name: 1-[(2S)-1-oxo-2-(3,4,5-trimethoxyphenyl)butyl]-2-piperidinecarboxylic acid, (2S,2'S)-2,2'-[1,2-ethanediylbis[imino(2-oxo-2,1-ethanediyl)oxy-3,1-phenylene[(1R)-3-(3,4-dimethoxyphenyl)propylidene]]] ester

Synonym: AP1903

MF: $C_{78}H_{98}N_4O_{20}$

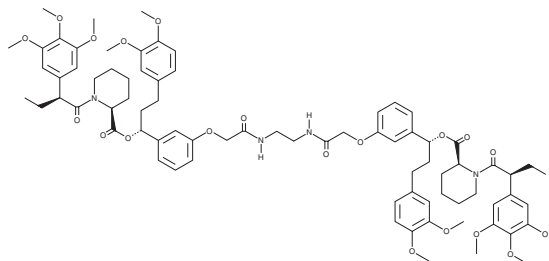
FW: 1,411.6

Purity: $\geq 90\%$ (mixture of isomers)

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

Rimiducid is supplied as a solid. A stock solution may be made by dissolving the rimiducid in the solvent of choice, which should be purged with an inert gas. Rimiducid is soluble in ethanol.

Description

Rimiducid is a chemical inducer of dimerization (CID) for fusion proteins of the FK506 binding protein (FKBP) mutant FKBP^{F36V}.¹ It selectively binds to FKBP^{F36V} over wild-type FKBP (IC_{50} s = 0.005 and $>10 \mu\text{M}$, respectively). Rimiducid induces apoptosis in HT-1080 fibrosarcoma cells expressing a FKBP^{F36V} fusion protein with the intracellular domain of Fas, also known as CD95, that contains a myristoylation sequence (EC_{50} = $\sim 0.1 \text{ nM}$). It also induces apoptosis in primary human T lymphocytes expressing a FKBP^{F36V}-Fas fusion protein (IC_{50} = $\sim 0.2 \text{ nM}$).² Rimiducid decreases serum human growth hormone (hGH) levels in mice injected with HT-1080 cells expressing FKBP^{F36V}-Fas that constitutively secrete hGH, indicating conditional cell ablation (ED_{50} = 0.4 mg/kg).¹

References

1. Clackson, T., Yang, W., Rozamus, L.W., *et al.* Redesigning an FKBP–ligand interface to generate chemical dimerizers with novel specificity. *Proc. Natl. Acad. Sci. USA* **95**(18), 10437-10442 (1998).
2. Thomis, D.C., Marktel, S., Bonini, C., *et al.* A Fas-based suicide switch in human T cells for the treatment of graft-versus-host disease. *Blood* **97**(5), 1249-1257 (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, 10/07/2024

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