

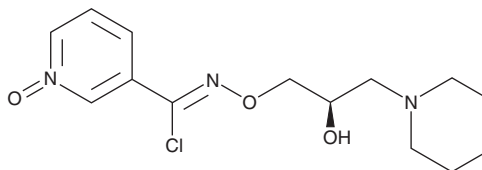
PRODUCT INFORMATION



Arimoclomol

Item No. 37368

CAS Registry No.: 289893-25-0
Formal Name: N-[(2R)-2-hydroxy-3-(1-piperidinyl)propoxy]-3-pyridinecarboximidoyl chloride, 1-oxide
Synonym: BRX-220
MF: $C_{14}H_{20}ClN_3O_3$
FW: 313.8
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Description

Arimoclomol is a co-inducer of heat shock proteins (Hsps).^{1,2} It increases pancreatic Hsp60 and Hsp70 levels, decreases malondialdehyde (MDA) and protein carbonyl levels, increases glutathione peroxidase (GPX) activity, and decreases Cu/Zn superoxide dismutase (SOD) activity in a rat model of cholecystokinin (CCK) octapeptide-induced acute pancreatitis when administered at a dose of 20 mg/kg.¹ Arimoclomol (10 mg/kg, i.p.) increases spinal cord levels of Hsp70 and Hsp90 and motoneuron survival, decreases disease severity, and increases lifespan in a Cu/Zn SOD mutant (SOD1^{G93A}) transgenic mouse model of amyotrophic lateral sclerosis (ALS).²

References

1. Rakonczay, Z., Jr., Iványi, B., Varga, I., *et al.* Nontoxic heat shock protein coinducer BRX-220 protects against acute pancreatitis in rats. *Free Radic. Biol. Med.* **32(12)**, 1283-1292 (2002).
2. Kieran, D., Kalmer, B., Dick, J.R.T., *et al.* Treatment with arimoclomol, a coinducer of heat shock proteins, delays disease progression in ALS mice. *Nat. Med.* **10(4)**, 402-405 (2004).

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, 12/21/2022

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