

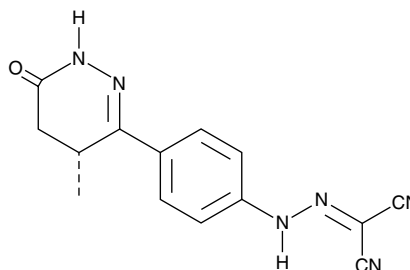
Product Information



Levosimendan

Item No. 16128

CAS Registry No.: 141505-33-1
Formal Name: 2-[2-[4-[(4R)-1,4,5,6-tetrahydro-4-methyl-6-oxo-3-pyridazinyl]phenyl]hydrazinylidene]-propanedinitrile
Synonyms: OR 1259, Simdax®, (R)-Simendan
MF: C₁₄H₁₂N₆O
FW: 280.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 250, 290, 382 nm



Laboratory Procedures

For long term storage, we suggest that levosimendan be stored as supplied at -20°C. It should be stable for at least two years.

Levosimendan is supplied as a crystalline solid. A stock solution may be made by dissolving the levosimendan in the solvent of choice. Levosimendan is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of levosimendan in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

Levosimendan is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Levosimendan is a calcium sensitizer that can cause increased cardiac contractility by binding troponin C (EC₅₀ = 9 nM), promotes vasodilation by activating ATP-sensitive potassium channels on vascular smooth muscle cells (EC₅₀ = 0.28 μM), and performs a cardioprotective function by prompting the opening of mitochondrial potassium channels in cardiomyocytes.¹⁻² It also has been reported to inhibit phosphodiesterases 3 and 4 in left ventricular cardiac tissue with IC₅₀ values of 2.5 nM and 25 μM, respectively.³

References

1. Nieminen, M.S., Fruhwald, S., Heunks, L.M.A., *et al.* Levosimendan: Current data, clinical use and future development. *Heart Lung Vessel* **5(4)**, 227-245 (2013).
2. Papp, Z., Csapó, K., Pollesello, P., *et al.* Pharmacological mechanisms contributing to the clinical efficacy of levosimendan. *Cardiovasc. Drug Rev.* **23(1)**, 71-98 (2005).
3. Szilgyi, S., Pollesello, P., Levijoki, J., *et al.* The effects of levosimendan and OR-1896 on isolated hearts, myocyte-sized preparations and phosphodiesterase enzymes of the guinea pig. *Eur. J. Pharmacol.* **486**, 67-74 (2004).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16128

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 07/11/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com