

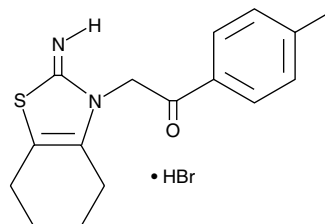
Product Information



Pifithrin- α

Item No. 13326

CAS Registry No.: 63208-82-2
Formal Name: 1-(4-methylphenyl)-2-(4,5,6,7-tetrahydro-2-imino-3(2H)-benzothiazolyl)-ethanone, monohydrobromide
Synonym: PFT- α
MF: C₁₆H₁₈N₂OS • HBr
FW: 367.3
Purity: $\geq 95\%$
Stability: ≥ 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 259 nm



Laboratory Procedures

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

Pifithrin- α is supplied as a crystalline solid. A stock solution may be made by dissolving the pifithrin- α in an organic solvent purged with an inert gas. Pifithrin- α is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pifithrin- α in these solvents is approximately 1 mg/ml.

Pifithrin- α is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pifithrin- α should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pifithrin- α has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that 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.

Pifithrin- α is an inactivator of p53 that blocks p53-dependent transcriptional activation and apoptosis.¹ It prevents p53-mediated apoptosis induced by cytotoxic compounds in C8 cells at $10 \mu\text{M}$ ¹ and in human umbilical vein endothelial cells at $30 \mu\text{M}$.² Pifithrin- α can also protect cells from DNA damage-induced apoptosis by a p53-independent mechanism that might involve cyclin D1.³

References

1. Komarov, P.G., Komarova, E.A., Kondratov, R.V., *et al.* A chemical inhibitor of p53 that protects mice from the side effects of cancer therapy. *Science* **285**, 1733-1737 (1999).
2. Lorenzo, E., Ruiz-Ruiz, C., Quesada, A.J., *et al.* Doxorubicin induces apoptosis and CD95 gene expression in human primary endothelial cells through a p53-dependent mechanism. *J. Biol. Chem.* **277**(17), 10883-10892 (2002).
3. Sohn, D., Graupner, V., Neise, D., *et al.* Pifithrin- α protects against DNA damage-induced apoptosis downstream of mitochondria independent of p53. *Cell Death and Differentiation* **16**, 869-878 (2009).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13326

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, 02/17/2015

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