

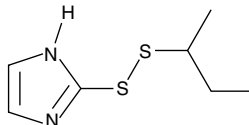
# Product Information



## PX 12

Item No. 14192

**CAS Registry No.:** 141400-58-0  
**Formal Name:** 2-[(1-methylpropyl)dithio]-1H-imidazole  
**MF:** C<sub>7</sub>H<sub>12</sub>N<sub>2</sub>S<sub>2</sub>  
**FW:** 188.3  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 235 nm



### Laboratory Procedures

For long term storage, we suggest that PX 12 be stored as supplied at -20°C. It should be stable for at least two years.

PX 12 is supplied as a crystalline solid. A stock solution may be made by dissolving the PX 12 in the solvent of choice. PX 12 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of PX 12 in these solvents is approximately 20, 14, and 16 mg/ml, respectively.

PX 12 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.

PX 12 is a competitive, irreversible inhibitor of thioredoxin 1 (Trx1), acting by binding to Cys<sup>73</sup> of this enzyme.<sup>1</sup> As Trx1 is overexpressed in certain cancers, PX 12 is effective in suppressing the growth of cancer cells, with growth inhibition in a panel of 60 human tumors significantly correlated with expression of Trx1 mRNA.<sup>2,3</sup> Through its effects on Trx1, PX 12 reduces hypoxia-induced HIF-1α protein levels (IC<sub>50</sub> = 7.2 μM), as well as expression of VEGF (IC<sub>50</sub> = 10.4 μM) and iNOS (IC<sub>50</sub> = 18.1 μM), culminating in attenuation of the proliferation of MCF-7 cells (IC<sub>50</sub> = 1.9 μM) and HT-29 cells (IC<sub>50</sub> = 2.9 μM) as well as reduced microvessel density in MCF-7 tumor xenografts.<sup>4</sup> PX 12 also blocks Trx-mediated activation of extracellular transglutaminase 2 (IC<sub>50</sub> = ~3 μM) and Trx-1-increased expression and translation of CYP1B1.<sup>5,6</sup>

### References

1. Oblong, J.E., Chantler, E.L., Gallegos, A., *et al.* Reversible inhibition of human thioredoxin reductase activity by cytotoxic alkyl 2-imidazolyl disulfide analogues. *Cancer Chemother. Pharmacol.* **34**, 434-438 (1994).
2. Powis, G. and Montfort, W.R. Properties and biological activities of thioredoxins. *Annu. Rev. Biophys. Biomol. Struct.* **30**, 421-455 (2001).
3. Mukherjee, A. and Martin, S.G. The thioredoxin system: A key target in tumour and endothelial cells. *Br. J. Radiol.* **81**, S57-S68 (2008).
4. Welsh, S.J., Williams, R.R., Birmingham, A., *et al.* The thioredoxin redox inhibitors 1-methylpropyl 2-imidazolyl disulfide and pleurotin inhibit hypoxia-induced factor 1α and vascular endothelial growth factor formation. *Mol. Cancer Ther.* **2(3)**, 235-243 (2003).
5. Jin, X., Stamnaes, J., Klöck, C., *et al.* Activation of extracellular transglutaminase 2 by thioredoxin. *J. Biol. Chem.* **286(43)**, 37866-37873 (2011).
6. Husbeck, B. and Powis, G. The redox protein thioredoxin-1 regulates the constitutive and inducible expression of the estrogen metabolizing cytochromes P450 1B1 and 1A1 in MCF-7 human breast cancer cells. *Carcinogenesis* **23(10)**, 1625-1630 (2002).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/14192](http://www.caymanchem.com/catalog/14192)

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

#### MATERIAL 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 Material 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, 04/25/2013

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)