

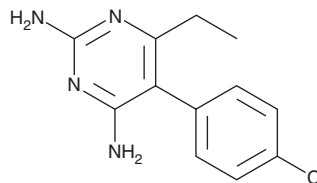
# PRODUCT INFORMATION



## Pyrimethamine

Item No. 16472

**CAS Registry No.:** 58-14-0  
**Formal Name:** 5,4-chlorophenyl-6-ethyl-2,4-pyrimidinediamine  
**Synonyms:** Darapram, Daraprim®, Khloridin, NSC 3061  
**MF:** C<sub>12</sub>H<sub>13</sub>ClN<sub>4</sub>  
**FW:** 248.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 285 nm  
**Supplied as:** A crystalline 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

Pyrimethamine is supplied as a crystalline solid. A stock solution may be made by dissolving the pyrimethamine in the solvent of choice. Pyrimethamine is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of pyrimethamine in these solvents is approximately 10 and 2.5 mg/ml, respectively.

Pyrimethamine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pyrimethamine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pyrimethamine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pyrimethamine is an antiprotozoal agent that is primarily active against *P. falciparum*, inhibiting the protozoal enzyme dihydrofolate reductase (DHFR).<sup>1</sup> By blocking DHFR activity, pyrimethamine prevents the production of tetrahydrofolic acid, an essential coenzyme involved in DNA and RNA synthesis. In *in vivo* antimalarial mouse models, pyrimethamine displays a prophylactic effect at an ED<sub>50</sub> value of 0.5 mg/kg.<sup>2</sup> Sulphonamides act synergistically with pyrimethamine by arresting production of dihydrofolic acid, resulting in the sequential blockade of the folate pathway of protozoa.<sup>3</sup> Pyrimethamine has also been found to limit the expression of the superoxide dismutase 1 (SOD1) gene, a metalloenzyme involved in amyotrophic lateral sclerosis.<sup>4</sup>

### References

1. Peters, W. Chemoprophylaxis and chemotherapy. *Brit. Med. J.* **2**, 95-98 (1971).
2. Adebajo, A.C., Odeiran, S.A., Aliyu, F.A., et al. *In vivo* antiplasmodial potentials of the combinations of four Nigerian antimalarial plants. *Molecules* **19**(9), 13136-13146 (2014).
3. Hwang, J., Bitarakwate, E., Pai, M., et al. Chloroquine or amodiaquine combined with sulfadoxine-pyrimethamine for uncomplicated malaria: A systematic review. *Trop. Med. Int. Health* **11**(6), 789-799 (2006).
4. Limpert, A.S., Mattmann, M.E., and Cosford, N.D.P. Recent progress in the discovery of small molecules for the treatment of amyotrophic lateral sclerosis (ALS). *Beilstein J. Org. Chem.* **9**, 717-732 (2013).

#### 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, 11/16/2022

#### 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