

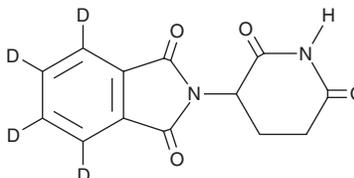
PRODUCT INFORMATION



(±)-Thalidomide-d₄

Item No. 29450

CAS Registry No.: 1219177-18-0
Formal Name: 2-(2,6-dioxo-3-piperidiny)-1H-isoindole-1,3(2H)-dione-4,5,6,7-d₄
Synonym: N-Phthaloylglutamide-d₄
MF: C₁₃H₆D₄N₂O₄
FW: 262.3
Chemical Purity: ≥98% ((±)-Thalidomide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A 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

(±)-Thalidomide-d₄ is intended for use as an internal standard for the quantification of (±)-thalidomide (Item No. 14610) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(±)-Thalidomide-d₄ is supplied as a solid. A stock solution may be made by dissolving the (±)-thalidomide-d₄ in the solvent of choice, which should be purged with an inert gas. (±)-Thalidomide-d₄ is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (±)-thalidomide-d₄ in these solvents is approximately 12 mg/ml.

Description

(±)-Thalidomide-d₄ is intended for use as an internal standard for the quantification of (±)-thalidomide (Item No. 14610) by GC- or LC-MS. (±)-Thalidomide is an immunomodulatory compound with diverse biological activities, including anticancer, anti-inflammatory, and teratogenic properties.¹⁻³ It prevents polymorphonuclear leukocyte (PMN) chemotaxis when used at concentrations of 1, 10, and 100 µg/ml.¹ (±)-Thalidomide increases IL-2-induced proliferation and IFN-γ production in primary human T cells *in vitro*.² It enhances natural killer (NK) cell-mediated cytotoxicity in MM.1S multiple myeloma cells.³ Thalidomide (4 mg/animal) reduces lung IL-6, TGF-β, VEGF, angiopoietin-1, angiopoietin-2, and collagen type Iα1 expression, inhibits pulmonary angiogenesis, and attenuates fibrosis in a mouse model of bleomycin-induced pulmonary fibrosis.⁴ It induces apoptosis in primary human embryonic fibroblasts (EC₅₀ = 8.9 µM) and induces limb and eye defects in chicken embryos (EC₅₀ = 50 µg/kg egg weight).⁵ Formulations containing thalidomide have been used in the treatment of multiple myeloma and erythema nodosum leprosum (ENL) in non-pregnant individuals.

References

1. Faure, M., Thivolet, J., and Gaucherand, M. *Arch. Dermatol. Res.* **269**(3), 275-280 (1980).
2. Haslett, P.A., Corral, L.G., Albert, M., *et al. J. Exp. Med.* **187**(11), 1885-1892 (1998).
3. Davies, F.E., Raje, N., Hideshima, T., *et al. Blood* **98**(1), 210-216 (2001).
4. Tabata, C., Tabata, R., Kadokawa, Y., *et al. J. Immunol.* **179**(1), 708-714 (2007).
5. Knobloch, J., Shaughnessy, J.D., Jr., and Rütther, U. *FASEB J.* **21**(7), 1410-1421 (2007).

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, 11/08/2022

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