

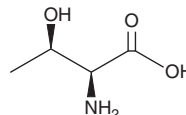
PRODUCT INFORMATION



L-Threonine

Item No. 30194

CAS Registry No.: 72-19-5
Synonym: (S)-Threonine
MF: $C_4H_9NO_3$
FW: 119.1
Purity: $\geq 95\%$
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

L-Threonine is supplied as a solid. Aqueous solutions of L-threonine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of L-threonine in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Threonine is an essential amino acid.¹ It undergoes catabolism via threonine dehydrogenase to yield the glucogenic metabolites succinyl coenzyme A (succinyl CoA) and pyruvate. Administration of L-threonine-deficient diets induces lameness of the thoracic and pelvic limbs and decreases food intake and weight gain in kittens.²

References

1. Massey, K.A., Blakeslee, C.H., and Pitkow, H.S. A review of physiological and metabolic effects of essential amino acids. *Amino Acids* **14**, 271-300 (1998).
2. Titchenal, C.A., Rogers, Q.R., Indrieri, R.J., *et al.* Threonine imbalance, deficiency and neurologic dysfunction in the kitten. *J. Nutr.* **110**(12), 2444-2459 (1980).

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, 10/26/2022

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