

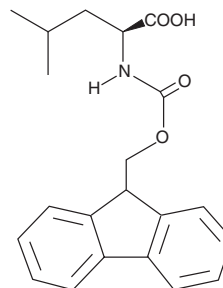
PRODUCT INFORMATION



Fmoc-L-Leucine

Item No. 10004888

CAS Registry No.: 35661-60-0
Formal Name: N-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-leucine
Synonyms: Fmoc-Leu, NPC 15199, NSC 334290
MF: C₂₁H₂₃NO₄
FW: 353.4
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 265, 289, 300 nm
Supplied as: A crystalline solid
Storage: Room Temperature
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Fmoc-L-leucine is supplied as a crystalline solid. A stock solution may be made by dissolving the Fmoc-L-leucine in the solvent of choice, which should be purged with an inert gas. Fmoc-L-leucine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Fmoc-L-leucine in these solvents is approximately 30 mg/ml.

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

Description

Peroxisome proliferator-activated receptor γ (PPAR γ) isoforms heterodimerize with retinoic X receptors to modulate gene expression related to adipocyte differentiation, fatty acid uptake and storage, and glucose metabolism.¹ Natural agonists of PPAR γ include fatty acids (e.g., linoleic acid and 15-deoxy- $\Delta^{12,14}$ -prostaglandin J₂), while thiazolidinediones (e.g., rosiglitazone and pioglitazone) are potent synthetic agonists.^{2,3} Fmoc-L-leucine is a partial agonist of PPAR γ .^{2,4} It activates PPAR γ with a lower potency ($K_i = 15$ versus 0.035 μ M) but a similar maximal efficacy compared to rosiglitazone.⁴ Fmoc-L-leucine improves insulin resistance in normal, diet-induced glucose-intolerant and in diabetic *db/db* mice, yet has reduced adipogenic activity.⁴ As a result, it is classified as a selective PPAR γ modulator (SPPARM), capable of producing insulin-sensitizing effects while minimizing side effects associated with full agonists.²

References

1. Heikkinen, S., Auwerx, J., and Argmann, C.A. PPAR γ in human and mouse physiology. *Biochim. Biophys. Acta* **1771**(8), 999-1013 (2007).
2. Villacorta, L., Schopfer, F.J., Zhang, J., et al. PPAR γ and its ligands: Therapeutic implications in cardiovascular disease. *Clin. Sci.* **116**(3), 205-218 (2009).
3. Zieleniak, A., Wójcik, M., and Wozniak, L.A. Structure and physiological functions of the human peroxisome proliferator-activated receptor γ . *Arch. Immunol. Ther. Exp.* **56**(5), 331-345 (2008).
4. Rocchi, S., Picard, F., Vamecq, J., et al. A unique PPAR γ ligand with potent insulin-sensitizing yet weak adipogenic activity. *Mol. Cell* **8**(4), 737-747 (2001).

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, 05/15/2023

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