

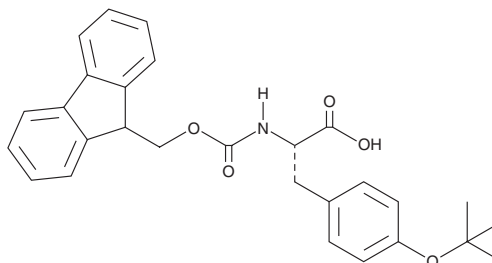
PRODUCT INFORMATION



Fmoc-Tyr(tBu)-OH

Item No. 42581

CAS Registry No.: 71989-38-3
Formal Name: O-(1,1-dimethylethyl)-N-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-tyrosine
Synonyms: Fmoc-O-tert-butyl-L-tyrosine, NSC 334301
MF: C₂₈H₂₉NO₅
FW: 459.5
Purity: ≥98%
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

Fmoc-Tyr(tBu)-OH is supplied as a solid. A stock solution may be made by dissolving the Fmoc-Tyr(tBu)-OH in the solvent of choice, which should be purged with an inert gas. Fmoc-Tyr(tBu)-OH is soluble (≥10 mg/ml) in DMSO and is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Fmoc-Tyr(tBu)-OH is an amino acid-containing building block.¹ It has commonly been used in solid-phase peptide synthesis (SPPS) to prevent tyrosine side-chain acylation.

Reference

1. Doherty-Kirby, A.L. and Lajoie, G.A. Solid-phase synthesis: A practical guide 129-196 (2000).

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, 03/13/2025

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