

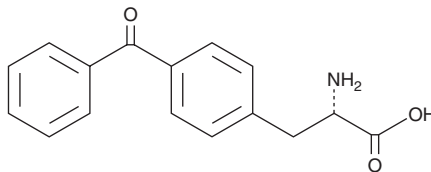
PRODUCT INFORMATION



p-Benzoyl-L-phenylalanine

Item No. 36168

CAS Registry No.: 104504-45-2
Formal Name: 4-benzoyl-L-phenylalanine
Synonym: BzF
MF: C₁₆H₁₅NO₃
FW: 269.3
Purity: ≥95%
UV/Vis.: λ_{max}: 259 nm
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

p-Benzoyl-L-phenylalanine (BzF) is supplied as a solid. A stock solution may be made by dissolving the BzF in the solvent of choice, which should be purged with an inert gas. BzF is soluble in hydrochloric acid.

Description

BzF is a photoreactive derivative of the amino acid L-phenylalanine (Item No. 31498).¹ It can be incorporated into synthetic peptides and used as a photolabel to cross-link the peptides to their protein binding partners to identify the residues involved in peptide-protein interactions. BzF has also been incorporated into G protein-coupled receptors (GPCRs) to identify potential ligand binding sites.²

References

1. Kauer, J.C., Erickson-Viitanen, S., Wolfe, H.R., Jr., et al. *p*-Benzoyl-L-phenylalanine, a new photoreactive amino acid. Photolabeling of calmodulin with a synthetic calmodulin-binding peptide. *J. Biol. Chem.* **261**(23), 10695-10700 (1986).
2. Becker, J.M. and Naider, F. Cross-linking strategies to study peptide ligand-receptor interactions. *Methods Enzymol.* **556**, 527-547 (2015).

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, 09/26/2022

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