

# PRODUCT INFORMATION



## Brain-Derived Basic Fibroblast Growth Factor (1-24) (bovine) (trifluoroacetate salt)

Item No. 24760

**Formal Name:** L-prolyl-L-alanyl-L-leucyl-L-prolyl-L- $\alpha$ -glutamyl-L- $\alpha$ -aspartylglycylglycyl-L-serylglycyl-L-alanyl-L-phenylalanyl-L-prolyl-L-prolylglycyl-L-histidyl-L-phenylalanyl-L-lysyl-L- $\alpha$ -aspartyl-L-prolyl-L-lysyl-L-arginyl-L-leucyl-L-tyrosine, trifluoroacetate salt

**Synonym:** Brain-Derived bFGF (1-24) (bovine)

**MF:**  $C_{118}H_{173}N_{31}O_{33} \cdot XCF_3COOH$

**FW:** 2,553.8

**Purity:**  $\geq 95\%$

**Supplied as:** A lyophilized powder

**Storage:**  $-20^\circ C$

**Stability:**  $\geq 4$  years

H-Pro—Ala—Leu—Pro—Glu—Asp—Gly—Gly—Ser—Gly—  
Ala—Phe—Pro—Pro—Gly—His—Phe—Lys—Asp—Pro—  
Lys—Arg—Leu—Tyr—OH  
 $\cdot XCF_3COOH$

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

### Laboratory Procedures

Brain-derived basic fibroblast growth factor (brain-derived bFGF) (1-24) (bovine) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the brain-derived bFGF (1-24) (bovine) (trifluoroacetate) in water. The solubility of brain-derived bFGF (1-24) (bovine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Brain-derived bFGF (1-24) is a peptide fragment of brain-derived bFGF.<sup>1</sup> bFGF is a peptide produced in bovine brain that is protective in a rat model of pressure-induced retinal ischemia. bFGF (1-24) corresponds to amino acid residues 1-24 of the full length peptide.

### Reference

1. Unoki, K. and LaVail, M.M. Protection of the rat retina from ischemic injury by brain-derived neurotrophic factor, ciliary neurotrophic factor, and basic fibroblast growth factor. *Invest. Ophthalmol. Vis. Sci.* **35**(3), 907-915 (1994).

#### 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, 11/17/2022

#### 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