

PRODUCT INFORMATION

Brain-Derived Acidic Fibroblast Growth Factor (102-111) (bovine) (trifluoroacetate salt)

Item No. 24761

Formal Name:	L-histidyl-L-alanyl-L- α -glutamyl-L-lysyl-L-histidyl-L-tryptophyl-L-phenylalanyl-L-valylglycyl-L-leucine, (trifluoroacetate salt)	H—His—Ala—Glu—Lys—His—Trp—Phe—Val—Gly—Leu—OH
Synonym:	aFGF (102-111)	
MF:	$C_{59}H_{82}N_{16}O_{13} \cdot XCF_3COOH$	$\cdot XCF_3COOH$
FW:	1,223.4	
Purity:	$\geq 95\%$	
Supplied as:	A lyophilized powder	
Storage:	-20°C	
Stability:	≥ 2 years	

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

Laboratory Procedures

Brain-derived acidic fibroblast growth factor (102-111) (bovine) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the brain-derived acidic fibroblast growth factor (102-111) (bovine) (trifluoroacetate salt) in water. The solubility of brain-derived acidic fibroblast growth factor (102-111) (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 acidic fibroblast growth factor (102-111) is a peptide fragment of brain-derived acidic fibroblast growth factor (aFGF).¹ aFGF is an angiogenic vascular endothelial cell mitogen produced in bovine brain that has sequence homology to interleukin-1.² It also shares sequence homology with the known neuropeptides neuromedin C (Item No. 24706), bombesin (Item No. 24151), neuromedin K (neurokinin B; Item No. 24542), substance K (neurokinin A; Item No. 24221), substance P (Item No. 24035), physalaemin, and eledoisin. aFGF (102-111) corresponds to amino acid residues 102-111 of the fulllength peptide.¹

References

1. Gimenez-Gallego, G., Rodkey, J., Bennett, C., *et al.* Brain-derived acidic fibroblast growth factor: Complete amino acid sequence and homologies. *Science* **230(4732)**, 1385-1388 (1985).
2. Thomas, K.A., Rios-Candelore, M., Giménez-Gallego, G., *et al.* Pure brain-derived acidic fibroblast growth factor is a potent angiogenic vascular endothelial cell mitogen with sequence homology to interleukin 1. *Proc. Natl. Acad. Sci. U.S.A.* **82(19)**, 6409-6413 (1985).

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