

# PRODUCT INFORMATION



## Bevonescein (trifluoroacetic acid)

Item No. 44659

**Formal Name:** N<sup>2</sup>-[(3',6'-dihydroxy-3-oxospiro[isobenzofuran-1(3H) 9'-[9H]xanthen]-5-yl)carbonyl]-L-glutaminy-L-valyl-L-prolyl-L-tryptophyl-L-α-glutamyl-L-α-glutamyl-L-prolyl-L-tyrosyl-L-tyrosyl-L-valyl-L-valyl-L-lysyl-L-lysyl-L-seryl-L-serylglycyl-glycinamide

**Synonym:** ALM-488

**Peptide Sequence:** 5-FAM-QVPWEEPYYVVKSSGG-NH<sub>2</sub>

**MF:** C<sub>112</sub>H<sub>144</sub>N<sub>22</sub>O<sub>32</sub> • XCF<sub>3</sub>COOH

**FW:** 2,310.5

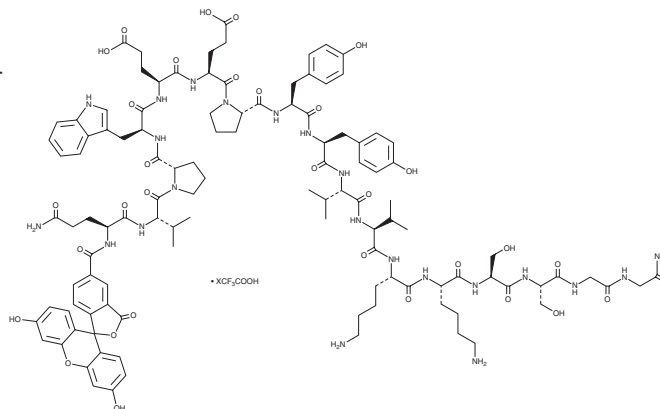
**Purity:** ≥95%

**Ex./Em. Max:** 488/530 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

Bevonescein (trifluoroacetic acid) is supplied as a solid. A stock solution may be made by dissolving the bevonescein (trifluoroacetic acid) in the solvent of choice, which should be purged with an inert gas. Bevonescein (trifluoroacetic acid) is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

### Description

Bevonescein is a peptide-dye conjugate for the visualization of intact and degenerated nerves.<sup>1</sup> It is composed of a synthetic 17-amino acid peptide conjugated to 5-carboxy fluorescein (5-FAM; Item No. 19581) and displays excitation/emission maxima of 488/530 nm.<sup>2,3</sup> Bevonescein (900 nmol/animal) has been used to identify degenerated nerves *via* real-time fluorescence imaging in mice in a model of facial nerve transection.<sup>1</sup>

### References

1. Crawford, K.L., Lee, Y.-J., Hom, M., *et al.* Identification of degenerated murine facial nerves with fluorescence labeling after transection injury. *Otolaryngol. Head Neck Surg.* **169(2)**, 234-242 (2023).
2. Lee, Y.J., Orosco, R.K., Bouvet, M., *et al.* Intraoperative nerve-specific fluorescence visualization in head and neck surgery: A Phase 1 trial. *Nat. Commun.* **16(1)**, 6060 (2025).
3. Azari, F., Zhang, K., Kennedy, G.T., *et al.* Precision surgery guided by intraoperative molecular imaging. *J. Nucl. Med.* **63(11)**, 1620-1627 (2022).

#### 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, 02/04/2026

#### 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