

# Product Information

## Biotin Cadaverine, Trifluoroacetate Salt

**Catalog Number:** 90063

**Unit Size:** 50 mg

### Storage and Handling

Store at -20°C or 4°C. Product is stable for at least 12 months from date of receipt when stored as recommended.

**Molecular Information:** C<sub>17</sub>H<sub>29</sub>F<sub>3</sub>N<sub>4</sub>O<sub>4</sub>S

**Molecular Weight:** 443

**Color and Form:** White solid

**Solubility:** Soluble in DMSO

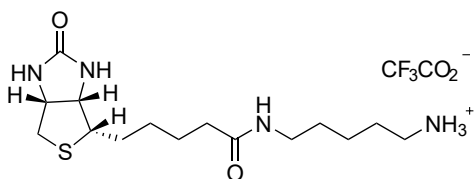


Figure 1. Biotin Cadaverine, Trifluoroacetate Salt

### Product Description

Biotin Cadaverine, Biotin-X Cadaverine (Cat. No. 90061), and Biotin Ethylenediamine (Cat. No. 90057) can be coupled to carboxylic acids, DNA and other biomolecules.

### Related Products

Cat. No.	Product
59002	EDC (EDAC)
90082	DMSO, Anhydrous
90050	Biotin SE
90051	Biotin-X, Succinimidyl Ester
90052	Biotin-XX, Succinimidyl Ester
90053	Biotin-X, Free Acid
90054	Biotin-XX, Free Acid
90057	Biotin Ethylenediamine, Hydrobromide (Neurobiotin™)
90058	Biotin-X-C5-Maleimide
90060	Biocytin Hydrazide
90061	Biotin-X Cadaverine, Trifluoroacetate Salt
90064 - 90066	Biotin MTSEA
90067	Biotin-PEO3-Amine
90068	Biotin-PEO4-Amine
90069	Biotin-PEO4-Propionate SE
90071	Biotin Ethylenediamine (free base)
90072	Biotin D(+)
90075	Biotin Ethylenediamine, Hydrochloride
90078	Biotin-PEO2-PPO2-Amine, TFA Salt
90079	Biotin Cadaverine, Free Base
90080	Biotin-X Cadaverine, Free Base
90083	Biotin-12 SE
90111	Biotin-PEO4-Hydrazide
90112	Biotin-PEO4-Maleimide
90113	Aminoxy-Biotin
92167	Biotin Azide
92168	Biotin Alkyne
92169	Biotin BCN
92186	Biotin Picolyl Azide
29086 - 29081	Streptavidin Conjugates
92035 - 96010	CF® Dye Amine
92208 - 92228	CF® Dye & Biotin SE Protein Labeling Kits

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.