



L-tyrosylglycyl-L-arginyl-L-lysyl-L-lysyl-L-arginyl-L-arginyl-L-glutaminyl-L-arginyl-L-arginyl-L-arginyl-L-lysyl-L-glutaminyl-L-isoleucyl-L- $\alpha$ -glutamyl-L-isoleucyl-L-lysyl-L-lysyl-L-phenylalanyl-L-lysine, trifluoroacetate salt

|                       |                           |                |                                          |
|-----------------------|---------------------------|----------------|------------------------------------------|
| Purity Specification: | ≥95%                      |                |                                          |
| Molecular formula:    | C119H212N46O26 • XCF3COOH |                |                                          |
| <b>CAS number:</b>    | <b>Formula weight:</b>    | <b>2703.25</b> | <b>Expiration date:</b> <b>09NOV2029</b> |

| Tests     | Results          |
|-----------|------------------|
| HPLC      | Purity: 100.0 %  |
| Mass Spec | [M-2H]2-: 1350.6 |

Gibby Perry

**WWW.CAYMANCHEM.COM**