



Item No. 24890 \* Batch No. 0526527

CAS number:                      Formula weight:    1513.68                      Expiration date:    10JUN2030

| Tests           | Results        |
|-----------------|----------------|
| Mass Spec       | M+: 1512.9     |
| HPLC            | Purity: 95.2 % |
| Peptide content | 0.00           |
| Peptide Co      | 81.00          |

Reviewed and approved by: Timothy Schrot

*Tim Sauer*

**WWW.CAYMANCHEM.COM**