

## FMOC-Statine [FMOC-Sta(3S,4S)-OH] \*CAS 158257-40-0\*

Catalog number: 5106, 5107 Unit size: 1 g, 10 g

| Product Details     |                          |
|---------------------|--------------------------|
| Storage Conditions  | Freeze (<-15 °C)         |
| Expiration Date     | 12 months upon receiving |
| Chemical Properties |                          |
| Appearance          | White solid              |
| Molecular Weight    | 397.47                   |
| Soluble In          | DMF                      |
| Chemical Structure  |                          |

## Applications

(3S,4S)-Statine is a novel amino acid of pepstatin, a low molecular weight inhibitor of acid proteases. This FMOC-Statine building lock is a very useful precursor for preparing protease-resistant peptide for pharmacetical application using the standard FMOC chemistry.