

## Product Information

**Description**

Biotin-PEG3-amine is reactive with NHS ester or carboxylic acid in the presence of EDC or HATU. PEG3 increases aqueous solubility of the molecules conjugated to the biotin compound. It also helps to minimize steric hindrance involved with the binding to avidin molecules.

**Molecular Formula:**

C<sub>18</sub>H<sub>34</sub>N<sub>4</sub>O<sub>5</sub>S

**Supplied As:**

Solid

**Stability:**

At least 2 years at -20°C. Avoid freeze/thaw cycles.

**Storage:**

-20°C

**MW:**

418.6 Da

**Cas No.**

359860-27-8

**Purity:**

≥98% by HPLC

## Chemical Structure

