

## Product Information

**Description**

tert-Butyl (6-iodohexyl)carbamate is a six carbon linker containing an iodine group and a t-butyl ester. The iodine (I) acts as a good leaving group for nucleophilic substitution reactions. The t-butyl ester groups can be deprotected under acidic conditions.

**Molecular Formula:**C<sub>11</sub>H<sub>22</sub>I NO<sub>2</sub>**Supplied As:**

Solid

**Stability:**

At least 6 months at -20°C. Avoid freeze/thaw cycles.

**Storage:**

-20°C

**MW:**

327.2 Da

**Cas No.**

172846-36-5

## Chemical Structure

