

Product Information

Description

5-Azidopentanoic acid ethyl ester is an azide compound that can be used to synthesize other compounds. 5-Azidopentanoic acid ethyl ester has been shown to inhibit the proliferation of cancer cells in vitro and in vivo. 5-Azidopentanoic acid ethyl ester reacts with amines to form a tethered compound, which can then be used as a probe for identifying amine-containing proteins.

Molecular Formula:C₇H₁₃N₃O₂**Supplied As:**

Solid

Stability:

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

Storage:

-20°C

MW:

171.2 Da

Cas No.

89896-39-9

Chemical Structure

