

Product Information

Description

Benzyl (5-hydroxypentyl)carbamate is a linker capable of repairing damaged DNA by binding to the 5'-hydroxyl group on the sugar backbone. It can also bind to DNA ligands and can inhibit cell proliferation in resistant cell lines. However, benzyl (5-hydroxypentyl)carbamate is toxic to cells and can also cause DNA damage and strand breakage.

Molecular Formula:

C₁₃H₁₉NO₃

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

237.3 Da

Cas No.

87905-98-4

Chemical Structure

