

Product Information

Description

MS4078 is a heterobifunctional linker that binds to protein targets and delivers cytotoxic drugs to cancer cells. It has a selective binding affinity for cancer cells and has shown minimal toxicity in pre-clinical studies.

Molecular Formula:

C₄₅H₅₂ClN₉O₈S

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

914.5 Da

Cas No.

2229036-62-6

Chemical Structure

