

Product Information

Description	MC-MMAF is a precursor of antibody drug conjugate. Mal group is reactive with thiol bearing molecule. MMAF is a synthetic antineoplastic agent.
Molecular Formula:	C ₄₉ H ₇₆ N ₆ O ₁₁
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	925.2 Da
Cas No.	863971-19-1
Purity:	98%

Chemical Structure

