

Product Information

Description	dBET1 is a PROTAC compound comprised of BET bromodomain antagonist (+)-JQ1 linked to a cereblon E3 ubiquitin ligase ligand.
Molecular Formula:	C ₃₈ H ₃₇ ClN ₈ O ₇ S
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	785.3 Da
Cas No.	1799711-21-9

Chemical Structure

