

## Product Information

<b>Description</b>	dBET6 is a potent, selective, and cell-permeable PROTAC connected by ligands for Cereblon and BET and has antitumor activity.
<b>Molecular Formula:</b>	C <sub>42</sub> H <sub>45</sub> ClN <sub>8</sub> O <sub>7</sub> S
<b>Supplied As:</b>	Solid
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	841.4 Da
<b>Cas No.</b>	1950634-92-0

## Chemical Structure

