

Product Information

Description

MS-436, through a set of water-mediated interactions, exhibits low nanomolar affinity (estimated K_i of 30-50 nM) with preference for the first bromodomain over the second. MS-436 effectively blocks BRD4 transcriptional activity in lipopolysaccharide-induced production of both nitric oxide and IL-6 in mouse macrophages (IC_{50} values are 3.8 and 4.9 μ M, respectively). MS-436 represents a new class of bromodomain inhibitors and will facilitate further investigation of the biological functions of the two bromodomains of BRD4 in gene expression.

Molecular Formula: $C_{18}H_{17}N_5O_3S$ **Supplied As:**

Solid

Stability:At least 6 months at $-20^{\circ}C$. Avoid freeze/thaw cycles.**Storage:** $-20^{\circ}C$ **MW:**

383.4 Da

Cas No.

1395084-25-9

Purity:

98%

Chemical Structure

