

(+)-JQ1

JQ1 is a chemical compound which is under development as a non-hormonal male contraceptive drug. Its chemical structure is a triazolothienodiazepine, a series originally discovered by Mitsubishi Pharma [WO/2009/084693], related to benzodiazepinederivatives such as etizolam, however JQ1 does not produce sedative or anxiolytic effects and is instead a potent and selective inhibitor of the bromodomain testis-specific protein BRDT, which is essential for chromatin remodeling during spermatogenesis.

Materials Provided

Catalog Number:	92-1149
Quantity Supplied:	1 vial(s), 1 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	457 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	1 mM
Reconstitution Volume:	2.188 mL
Reconstitution Solvent:	DMSO

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com