

## RVX-208

**Catalog #:** 27111

**Lot #:** 140113

**Size:** 5 mg

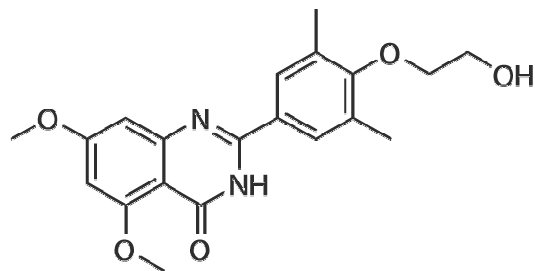
**Structure:**

**CAS Registry #:** 1044870-39-4

**Purity:** ≥98%

**Chemical Formula:** C<sub>20</sub>H<sub>22</sub>N<sub>2</sub>O<sub>5</sub>

**Molecular Weight:** 370.4



**Description:** RVX-028, also known as RVX-000222, is a potent BET bromodomain inhibitor. It preferentially binds to the second bromodomain on BET proteins. It also stimulates apolipoprotein (APO) AI gene expression, which increases apoA-I and HDL-C *in vitro*. *In vivo*, RVX-208 enhances cholesterol efflux via different pathways and significantly increases serum apoA-I and HDL-C in monkeys.

**Appearance:** A crystalline solid

**Solubility:** Solubility in DMSO is 74 mg/ml. Solubility in water and ethanol is <1 mg/ml.

**Biological Activity:** RVX-208 has been shown to have an IC<sub>50</sub> = 0.510 μM for BD2, which is ~170-fold selectivity over BD1.

**Storage/Stability:** Store at or below -20°C for up to two years.

**Quality Control:** The purity was determined by HPLC analysis.

### References:

1. Bailey, D., *et al.*, *J Am Coll Cardiol.* 2010;**55(23)**:2580-2589.
2. Picaud, S., *et al.*, *Proc Natl Acad Sci U.S.A.* 2013; **110(49)**:19754-19759.
3. McNeill, E. *Curr Opin Investig Drugs.* 2010;**11(3)**:357-364.