

RVX-208

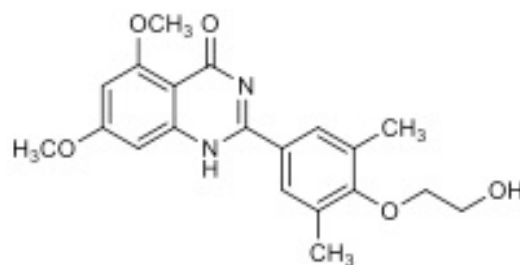
Catalog No: 14090, 14091**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 370.4

C₂₀H₂₂N₂O₅

CAS 1044870-39-4

Physical Properties: White powder

Names: RVX-208, 2-[4-(2-hydroxyethoxy)-3,5-dimethylphenyl]-5,7-dimethoxy-4(3H)-quinazolinone

Pharmacology: RVX-208 is a potent BET bromodomain antagonist with (IC₅₀ = 0.51 µM for BD2) with approximately 170-fold selectivity over BD1 (ref 1,2). Increases apolipoprotein A-1 and HDL cholesterol *in vitro* and *in vivo* (ref 3). Reduces atherosclerosis (ref 4).

Solubilization: May be dissolved in DMSO (75 mg/ml); Ethanol (4 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. S Picaud *et al. Proc. Natl. Acad. Sci. USA* 2013, 110:19754
2. KG McLure *et al. PLoS One* 2013, 8:e83190
3. D Bailey *et al. J. Am. Coll. Cardiol.* 2010, 55:2580
4. R Jahagirdar *et al. Atherosclerosis* 2014, 236:91

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.