

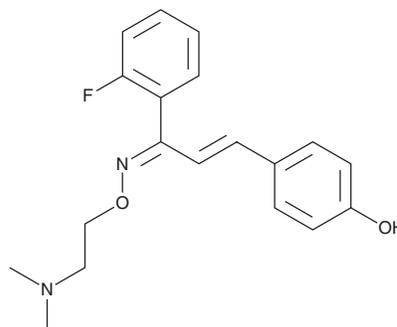
# PRODUCT INFORMATION



## Eplivanserin Mixture

Item No. 43749

**CAS Registry No.:** 130581-13-4  
**Formal Name:** 1-(2E-fluorophenyl)-3-(4-hydroxyphenyl)-2-propen-1-one, O-[2-(dimethylamino)ethyl]oxime  
**Synonyms:** SR 46349/SR 46615, SR 46349 and SR 46615 Mixture, SR 46349B/SR 46615, SR 46349B and SR 46615 Mixture  
**MF:** C<sub>19</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>2</sub>  
**FW:** 328.4  
**Purity:** ≥98% (mixture)  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Eplivanserin mixture is supplied as a solid. A stock solution may be made by dissolving the eplivanserin mixture in the solvent of choice, which should be purged with an inert gas. Eplivanserin mixture is soluble in organic solvents such as ethanol and DMSO. Eplivanserin mixture is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

### Description

Eplivanserin mixture is a mixture of the serotonin (5-HT) receptor subtype 5-HT<sub>2A</sub> antagonist SR 46349 and its isomer SR 46615.<sup>1,2</sup>

### References

1. Rinaldi-Carmona, M., Congy, C., Santucci, V., *et al.* Biochemical and pharmacological properties of SR 46349B, a new potent and selective 5-hydroxytryptamine<sub>2</sub> receptor antagonist. *J. Pharmacol. Exp. Ther.* **262**(2), 759-768 (1992).
2. Yang, L.H., Callas, D., Herbert, J.M., *et al.* Modulation of platelet function and vascular smooth muscle contractile actions by a novel, selective, highly potent 5-HT<sub>2</sub> antagonist (SR46349). *Thromb. Res.* **81**(5), 555-567 (1996).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/16/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM