

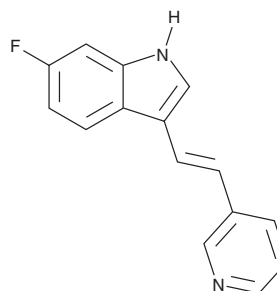
# PRODUCT INFORMATION



**680C91**

Item No. 34823

**CAS Registry No.:** 163239-22-3  
**Formal Name:** 6-fluoro-3-[(1E)-2-(3-pyridinyl)ethenyl]-1H-indole  
**MF:** C<sub>15</sub>H<sub>11</sub>FN<sub>2</sub>  
**FW:** 238.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 340 nm  
**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

680C91 is supplied as a solid. A stock solution may be made by dissolving the 680C91 in the solvent of choice, which should be purged with an inert gas. 680C91 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 680C91 in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

680C91 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 680C91 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 680C91 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

680C91 is an inhibitor of tryptophan 2,3-dioxygenase (TDO; IC<sub>50</sub>s = 1.46 and 0.28 μM for the mouse and human enzymes, respectively).<sup>1</sup> It is selective for TDO over indoleamine 2,3-dioxygenase (IDO; IC<sub>50</sub> = >80 μM for the mouse enzyme). 680C91 (40 μM) inhibits the migration of SK-MEL-28 melanoma cells induced by dexamethasone (Item No. 11015).<sup>2</sup> It increases brain levels of tryptophan, serotonin (5-HT), and 5-hydroxy indole-3-acetic acid (5-HIAA) in rats in a dose-dependent manner.<sup>3</sup>

## References

1. Pilotte, L., Larrieu, P., Stroobant, V., *et al.* Reversal of tumoral immune resistance by inhibition of tryptophan 2,3-dioxygenase. *Proc. Natl. Acad. Sci. USA* **109**(7), 2497-2502 (2012).
2. Cecchi, M., Paccosi, S., Silvano, A., *et al.* Dexamethasone induces the expression and function of tryptophan-2-3-dioxygenase in SK-MEL-28 melanoma cells. *Pharmaceuticals (Basel)* **14**(3), 211 (2021).
3. Salter, M., Hazelwood, R., Pogson, C.I., *et al.* The effects of a novel and selective inhibitor of tryptophan 2,3-dioxygenase on tryptophan and serotonin metabolism in the rat. *Biochem. Pharmacol.* **49**(10), 1435-1442 (1995).

### 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, 01/03/2023

## 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](http://WWW.CAYMANCHEM.COM)