

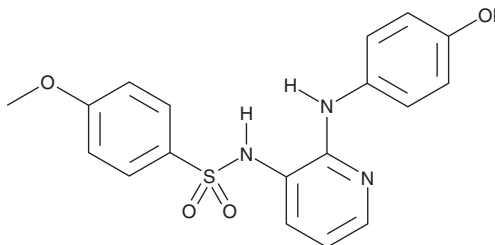
# PRODUCT INFORMATION



## ABT-751

Item No. 22935

**CAS Registry No.:** 141430-65-1  
**Formal Name:** N-[2-[(4-hydroxyphenyl)amino]-3-pyridinyl]-4-methoxybenzenesulfonamide  
**MF:** C<sub>18</sub>H<sub>17</sub>N<sub>3</sub>O<sub>4</sub>S  
**FW:** 371.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 244, 280 nm  
**Supplied as:** A crystalline 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

ABT-751 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABT-751 in the solvent of choice, which should be purged with an inert gas. ABT-751 is soluble in DMSO.

### Description

ABT-751 is a tubulin polymerization inhibitor that binds to the colchicine binding site on β-tubulin (K<sub>i</sub> = 3.3 μM).<sup>1</sup> It inhibits tubulin polymerization in a concentration-dependent manner and inhibits growth of 26 human tumor cell lines (IC<sub>50</sub>s = 0.06-0.08 μg/ml). ABT-751 (100 mg/kg, p.o.) induces tumor regression in NB-1382 neuroblastoma, IRS56 rhabdomyosarcoma, and KT-6 Wilms mouse xenograft models.<sup>2</sup> It also induces tumor endothelial cell retraction and reduces tumor perfusion, but not muscle perfusion, in a 9L rat glioma subcutaneous tumor model.<sup>3</sup>

### References

1. Yoshimatsu, K., Yamaguchi, A., Yoshino, H., *et al.* Mechanism of action of E7010, an orally active sulfonamide antitumor agent: Inhibition of mitosis by binding to the colchicine site of tubulin. *Cancer Res.* **57(15)**, 3208-3213 (1997).
2. Morton, C.L., Favours, E.G., Mercer, K.S., *et al.* Evaluation of ABT-751 against childhood cancer models in vivo. *Invest. New. Drugs* **25(4)**, 285-295 (2007).
3. Luo, Y., Hradil, V.P., Frost, D.J., *et al.* ABT-751, a novel tubulin-binding agent, decreases tumor perfusion and disrupts tumor vasculature. *Anticancer Drugs* **20(6)**, 483-492 (2009).

#### 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, 10/25/2022

#### 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