

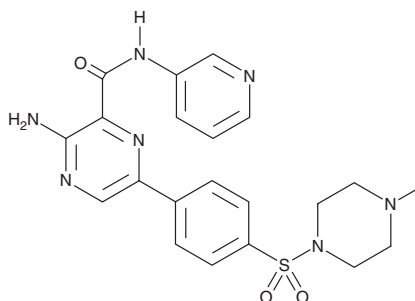
# PRODUCT INFORMATION



## AZD 2858

Item No. 16728

**CAS Registry No.:** 486424-20-8  
**Formal Name:** 3-amino-6-[4-[[4-methyl-1-piperazinyl)sulfonyl]phenyl]-N-3-pyridinyl-2-pyrazinecarboxamide  
**MF:** C<sub>21</sub>H<sub>23</sub>N<sub>7</sub>O<sub>3</sub>S  
**FW:** 453.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 304, 383 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

AZD 2858 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 2858 in the solvent of choice, which should be purged with an inert gas. AZD 2858 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of AZD 2858 in these solvents is approximately 5 and 10 mg/ml, respectively.

AZD 2858 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 2858 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AZD 2858 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Glycogen synthase kinase 3β (GSK3β) was originally identified for its ability to constitutively phosphorylate and inactivate glycogen synthase, preventing glycogen synthesis.<sup>1</sup> It can also phosphorylate proteins that are relevant to osteogenesis and Alzheimer's disease, the latter for which it has earned the moniker, tau phosphorylating kinase.<sup>2-4</sup> AZD 2858 is a pyrazine analog that inhibits GSK3β with a K<sub>i</sub> value of 4.9 nM.<sup>5</sup> It crosses the blood brain barrier and inhibits tau phosphorylation *in vitro* with an IC<sub>50</sub> value of 76 nM.<sup>5</sup> AZD 2858 has also been shown to increase bone mass (*via* Wnt activation) in rats after a two-week treatment with a maximum effective oral dose of 20 mg/kg once daily.<sup>6</sup> Furthermore, by inhibiting GSK3β, AZD 2858 can stabilize β-catenin in rat mesenchymal stem cells (EC<sub>50</sub> = 234 nM), spurring osteoblast differentiation.<sup>7</sup>

### References

1. Ring, D.B., Johnson, K.W., Henriksen, E.J., *et al. Diabetes* **52**(3), 588-95 (2003).
2. Asuni, A.A., Hooper, C., Reynolds, C.H., *et al. Eur. J. Neurosci.* **24**(12), 3387-3392 (2006).
3. Phiel, C.J., Wilson, C.A., Lee, V.M., *et al. Nature* **423**(6938), 435-439 (2003).
4. Guerrero, F., Herencia, C., Almadén, Y., *et al. PLoS One* **9**(2), e89179 (2014).
5. Berg, S., Bergh, M., Hellberg, S., *et al. J. Med. Chem.* **55**(21), 9107-9119 (2012).
6. Marsell, R., Sisask, G., Nilsson, Y., *et al. Bone* **50**(3), 619-627 (2012).
7. Gilmour, P.S., O'Shea, P.J., Fagura, M., *et al. Toxicol. Appl. Pharmacol.* **272**(2), 399-407 (2013).

#### 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, 11/30/2022

#### 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