

# PRODUCT INFORMATION



## WDR77 (human, recombinant)

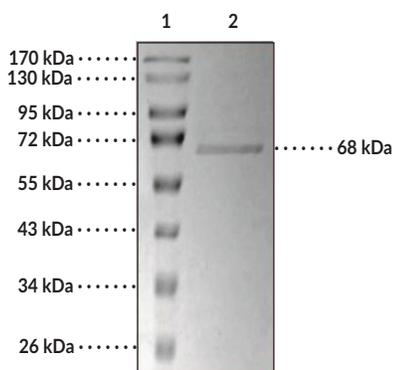
Item No. 43756

### Overview and Properties

<b>Synonyms:</b>	Androgen Receptor Cofactor p44, MEP-50, Methylosome Protein 50, Methylosome Protein WDR77, WD Repeat-containing protein 77
<b>Source:</b>	Recombinant human N-terminal GST-tagged WDR77 expressed in insect cells
<b>Amino Acids:</b>	1-342
<b>Uniprot No.:</b>	Q9BQA1
<b>Molecular Weight:</b>	68 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥90% estimated by SDS-PAGE
<b>Supplied in:</b>	Sterile 50 mM Tris-HCl, pH 7.5, with 50-300 mM sodium chloride, 10 mM glutathione, 0.1 mM EDTA, 0.25-1.0 mM DTT, 0-0.1 mM PMSF, and 10-25% glycerol
<b>Bioactivity:</b>	See figures for details

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

### Image



Lane 1: MW Markers  
Lane 2: WDR77

SDS-PAGE Analysis of WDR77. This protein has a calculated molecular weight of 68 kDa.

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, 09/25/2025

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

# PRODUCT INFORMATION



## Description

---

Methylosome protein WDR77, also known as methylosome protein 50 (MEP50) and androgen receptor cofactor p44, is an obligate cofactor of protein arginine methyltransferase 5 (PRMT5) and androgen receptor coactivator.<sup>1,2</sup> It contains 342 amino acid residues and six putative WD-40 repeats that act as scaffolds for protein-protein interactions.<sup>3</sup> WDR77 is ubiquitously expressed and localizes to both the cytoplasm and nucleus.<sup>1,2</sup> It binds to PRMT5 to stabilize and increase its activity, as well as recruits substrates to the PRMT5 complex for methylation.<sup>1</sup> WDR77 acts as a coactivator of androgen receptor transcription mediated through an androgen receptor-Smad1-WDR77 complex.<sup>2</sup> High cytoplasmic and low nuclear levels of WDR77 are associated with the development of castration-resistant prostate cancer. Cayman's WDR77 (human, recombinant) protein has a calculated molecular weight of 68 kDa.

## References

---

1. Antonyamy, S. The structure and function of the PRMT5:MEP50 complex. *Macromolecular protein complexes* 185-194 (2017).
2. Daniels, G., Jha, R., Shen, Y., *et al.* Androgen receptor coactivators that inhibit prostate cancer growth. *Am. J. Clin. Exp. Urol.* **5(1)**, 62-70 (2014).
3. Abe, Y., Suzuki, Y., Kawamura, K., *et al.* MEP50/PRMT5-mediated methylation activates GLI1 in Hedgehog signalling through inhibition of ubiquitination by the ITCH/NUMB complex. *Commun. Biol.* **2**, 23 (2019).

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