

# PRODUCT INFORMATION

## MR1/ $\beta$ 2-Microglobulin Complex (human, recombinant)

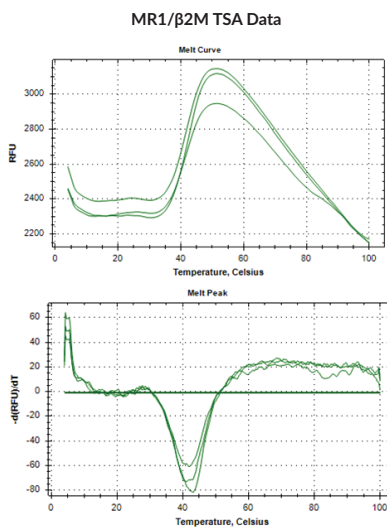
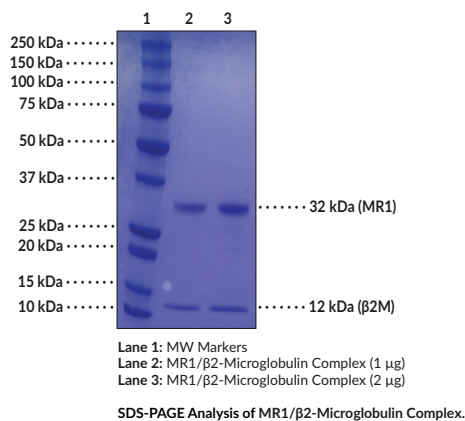
Item No. 30585

### Overview and Properties

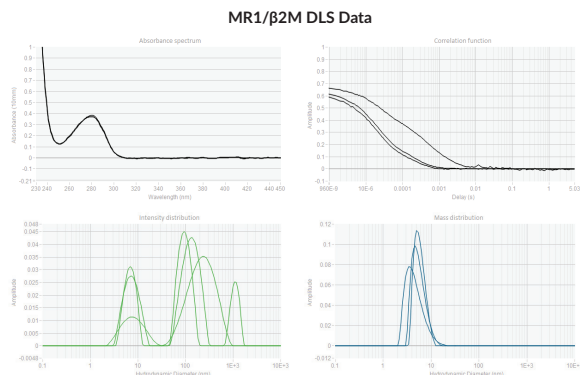
<b>Synonyms:</b>	<b>MR1:</b> Class I Histocompatibility Antigen-like Protein, Major Histocompatibility Complex Class I-related Gene Protein, MHC Class I-related Gene Protein <b><math>\beta</math>2-Microprotein:</b> $\beta$ 2M
<b>Source:</b>	Recombinant human MR1 and $\beta$ 2-microglobulin expressed in <i>E. coli</i>
<b>Amino Acids:</b>	22-292 (MR1), 21-119 ( $\beta$ 2-microglobulin)
<b>Uniprot No.:</b>	Q95460 (MR1), P61769 ( $\beta$ 2-microglobulin)
<b>Molecular Weight:</b>	32 (MR1) and 12 ( $\beta$ 2-microglobulin) kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	$\geq 1$ year
<b>Purity:</b>	$\geq 90\%$ estimated by SDS-PAGE
<b>Supplied in:</b>	50 mM HEPES, pH 8.0, with 150 mM sodium chloride
<b>Protein</b>	
<b>Concentration:</b>	<i>batch specific</i> mg/ml

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

### Images



Thermal shift assay (TSA) shows that the MR1/ $\beta$ 2M is well folded with a  $T_m$  of 43°C. These are representative data.



Dynamic light scattering (DLS) analysis showed that 99.8% of the protein by mass formed a complex with an average diameter of 27 nm. The Pdl of this complex was 0.32. These are representative data.

**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, 04/13/2023

**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

# PRODUCT INFORMATION



## Description

MHC class I-related gene protein (MR1) is a non-polymorphic MHC class Ib antigen-presenting cell surface molecule that is required for T cell receptor-mediated activation of mucosal-associated invariant T (MAIT) cells.<sup>1,2</sup> MR1 is composed of  $\alpha 1$  and  $\alpha 2$  domains, which form an antigen-binding pocket, and an  $\alpha 3$  domain that interacts with  $\beta 2$ -microglobulin, the light chain component of MHC class I molecules that facilitates complex transport to the cell surface and antigen presentation to cytotoxic T cells.<sup>2,3</sup> Upon binding of a microbial antigen, MR1 undergoes a conformational change in the endoplasmic reticulum and translocates to the cell surface with  $\beta 2$ -microglobulin where it induces MAIT cell activation *via* an interaction with the MAIT cell T cell receptor and activates various immunomodulatory effects, including cytokine release, initiation of adaptive immune responses, and promotion of tissue repair.<sup>4,5</sup> The antigen-binding domain of MR1 binds to vitamin B metabolites, including metabolites of vitamin B<sub>2</sub>, also known as riboflavin, and vitamin B<sub>9</sub> (folic acid; Item No. 20515).<sup>6</sup> MHC class I-associated  $\beta 2$ -microglobulin exhibits equilibrium exchange with circulating soluble  $\beta 2$ -microglobulin and serum levels of  $\beta 2$ -microglobulin are increased in patients with cancer, renal failure, systemic amyloidosis, or various autoimmune diseases, including multiple sclerosis.<sup>3,7-9</sup>

## References

1. Lamichhane, R. and Ussher, J.E. Expression and trafficking of MR1. *Immunology* **151**(3), 270-279 (2017).
2. Krovi, S.H. and Gapin, L. Structure and function of the non-classical major histocompatibility complex molecule MR1. *Immunogenetics* **68**(8), 549-559 (2016).
3. Nomura, T., Huang, W.-C., Zhau, H.E., *et al.*  $\beta 2$ -Microglobulin-mediated signaling as a target for cancer therapy. *Anticancer Agents Med. Chem.* **14**(3), 343-352 (2014).
4. McWilliam, H.E.G., Eckle, S.B.G., Theodossis, A., *et al.* The intracellular pathway for the presentation of vitamin B-related antigens by the antigen-presenting molecule MR1. *Nat. Immunol.* **17**(5), 531-537 (2016).
5. McWilliam, H.E.G. and Salio, M. Understanding and modulating the MR1 metabolite antigen presentation pathway. *Mol. Immunol.* **129**, 121-126 (2021).
6. Kjer-Nielsen, L., Patel, O., Corbett, A.J., *et al.* MR1 presents microbial vitamin B metabolites to MAIT cells. *Nature* **491**(7426), 717-723 (2012).
7. Eichner, T. and Radford, S.E. Understanding the complex mechanisms of  $\beta 2$ -microglobulin amyloid assembly. *FEBS J.* **278**(20), 3868-3883 (2011).
8. Stoppini, M. and Bellotti, V. Systemic amyloidosis: Lessons from  $\beta 2$ -microglobulin. *J. Biol. Chem.* **290**(16), 9951-9958 (2015).
9. Bagnato, F., Durastanti, V., Finamore, L., *et al.* Beta-2 microglobulin and neopterin as markers of disease activity in multiple sclerosis. *Neurol. Sci.* **24**(Suppl 5), S301-S304 (2003).

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