



Product Data Sheet

Product Name: IRBP, Interphotoreceptor Retinoid Binding Protein Fragment

Catalog Number: AS-60183-1 (1 mg) Lot Number: See label on vial
AS-60183-5 (5 mg)

Sequence: H-Ser-Gly-Ile-Pro-Tyr-Ile-Ile-Ser-Tyr-Leu-His-Pro-Gly-Asn-Thr-Ile-Leu-His-Val-Asp-OH (3-letter code)
SGIPYIISYLHPGNTILHVD (1-letter code)

Molecular Weight: 2210.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 µl to 1 mg IRBP peptide.

Storage: IRBP peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: This 20-residue peptide, a major pathogenic T-cell epitope (161–180), is present in the first homologous repeat of the interphotoreceptor retinoid binding protein peptide (IRBP). It has been shown to induce posterior uveitis (EAU). Ref: Sonoda, K. et al. *J. Immunol.* **171**, 2652 (2003); Read, RW. and SR. Barnum, *ARVO Meeting Abstracts* **43**, 2281 (2002); Silver, P. et al. *Invest. Ophthalmol. Vis. Sci.* **36**, 946 (1995); Karabekian, Z. et al. *Invest. Ophthalmol. Vis. Sci.* **46**, 3769 (2005).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

The Ag, a 20-aa peptide sequence SGIPYIISYLHPGNTILHVD encoded by human interphotoreceptor retinoid binding protein (IRBP) residues 161–180 (p161–180), encoding the major pathogenic epitope, was synthesized by AnaSpec using Fmoc technology. Ag pulsing was for 4.5 h, unless otherwise specified. For maturation the medium was additionally supplemented with 100 µg/ml p161–180 and/or 1 µg/ml LPS (*E. coli* O111:B4; Sigma-Aldrich), 5 µg/ml anti-CD40 (3/23; BD Biosciences)- [Tan, Jung. et al. *The Journal of Immunology.* **178**, 5578 \(2007\).](#)

Published Citations:

Tan, Jung. et al. *The Journal of Immunology.* **178**, 5578 (2007).

Related Product:

Name	Cat #	Size
IRBP (1-20), human (GPTHLFQPSLVLDMAKVLLD)	AS-62297	1 mg

For Research Use Only