

### HLA-B\*39:06 easYmer kit

<b>Catalog#</b>	1080-01
<b>Size</b>	20, 50, 150, or 500 tests
<b>easYmer</b>	HLA-B*39:06 (C67S; unpaired Cysteine in position 67 is substituted for Serine) - peptide receptive, biotinylated in Tris/Maleate pH 7 with 30% Glycerol
<b>Folding buffer</b>	Tris/Maleate pH 7
<b>Peptide</b>	RRYTRRISL, a stable HLA-B*39:06 binder Positive control for evaluation analysis of peptide-HLA folding
<b>Application</b>	easYmers can be used to generate monomers with your choice of peptide. The monomers can easily be tetramerized with fluorophore conjugated streptavidin and used to analyse T cells by flow cytometry. The easYmer reagent can also be used to evaluate specific peptide-HLA I binding.

**For Research Use Only (RUO)**

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) info@eaglebio.com



**EAGLE**  
BIOSCIENCES