

HLA-A*68:02 easYmer kit

Catalog#	1043-01
Size	20, 50, 150, or 500 tests
easYmer	HLA-A*68:02 - peptide receptive, biotinylated in Tris/Maleate pH 7 with 30% Glycerol
Folding buffer	Tris/Maleate pH 7
Peptide	ETVAIDRPA, a stable HLA-A*68:02 binder Positive control for evaluation analysis of peptide-HLA folding
Application	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 info@eaglebio.com



EAGLE
BIOSCIENCES