

HLA-C*07:02 (C1G) easYmer kit

| | |
|-----------------------|---|
| Catalog# | 1122-01 |
| Size | 20, 50, 150, or 500 tests |
| easYmer | HLA-C*07:02 (C1G; unpaired Cysteine in position 1 is substituted for Glycine) - peptide receptive, biotinylated in Tris/Maleate pH 7 with 30% Glycerol |
| Folding buffer | Tris/Maleate pH 7 |
| Peptide | IE1 309-317 (CRVLCCYVL), a known HLA-C*07:02 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)