

hERG Cardiac Safety Assay Service

Biochemical hERG Assay to assess cardiac safety of new compounds using fluorescence polarization for measuring compound binding to hERG.

CAT#:	hERG								
Target Overview									
Enzyme	Human ERG								
Probe:	Predictor Tracer, 1 nM (ThermoFisher)								
Assay Properties									
Readout	Competition binding assay with Tracer								
Reference Compound IC50s:	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>E-4031</td> <td>7.2</td> </tr> <tr> <td>Astemizole</td> <td>0.84</td> </tr> <tr> <td>GBR 12909</td> <td>21.5</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	E-4031	7.2	Astemizole	0.84	GBR 12909	21.5
Compounds	IC50 (nM)								
E-4031	7.2								
Astemizole	0.84								
GBR 12909	21.5								
Scientific Information:	P07900 (http://www.uniprot.org/uniprot/P07900)								

Control Compound IC50 for hERG

