

hERG

(Potassium voltage-gated channel subfamily H member 2, Ether-a-go-go-related gene potassium channel 1, ERG, ERG1, HERG, KCNH2)

- **CAT#:** hERG
- **Enzyme:** Human ERG
- **Probe:** Predictor™ Tracer, 1 nM
- **Reaction:**
Competition binding assay with Tracer.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
E-4031	7.2
Astemizole	0.84
GBR 12909	21.5

- **Scientific Information:** P07900 (<http://www.expasy.org/uniprot/P07900>)

Control Compound IC50 for hERG

