

---

## ARG(ABL2)

Product code: 08-102

Full-length human ARG [2-1161(end) amino acids of accession number NP\_001161708.1] was expressed as N-terminal GST-fusion protein (153 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : ABLtide

ATP ( $\mu$ M) Km app / Bin : 24 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 27

IC50 at 1 mM ATP (nM) : 400

---