

---

---

## EGFR[D770\_N771insNPG]

Product code: 08-553

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and D770\_N771insNPG of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system.

|                             |                 |
|-----------------------------|-----------------|
| Assay platform              | : MSA           |
| Substrate                   | : Srctide       |
| ATP ( $\mu$ M) Km app / Bin | : 2.3 / 5       |
| Metal                       | : Mg+Mn         |
| Reference compound          | : Staurosporine |
| IC50 at ATP Bin (nM)        | : 18            |
| IC50 at 1 mM ATP (nM)       | : 930           |