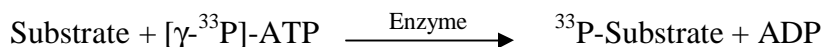


RAF1

(RAF proto-oncogene serine/threonine-protein kinase, Raf-1, C-RAF, cRaf, RAF)

- **CAT#:** RAF1
- **Enzyme:** Human RAF1
- **Substrate:** MEK1 (K97R), 0.5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|-----------|-----------|
| GW 5074 | 14 |

- **Scientific Information:** P04049 (<http://www.expasy.org/uniprot/P04049>)

Reference Compound IC50 for RAF1

