

ARAF

(B-Raf proto-oncogene serine/threonine-protein kinase, A-raf-1, Proto-oncogene Pks, ARAF1, PKS, PKS2)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
GW 5074	11.6

- **Scientific Information:** P10398 (<http://www.expasy.org/uniprot/P10398>)

Reference compound IC50 for ARAF

