

# EGFR (D770GY) Kinase Assay Service

Activity-based EGFR (D770GY) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

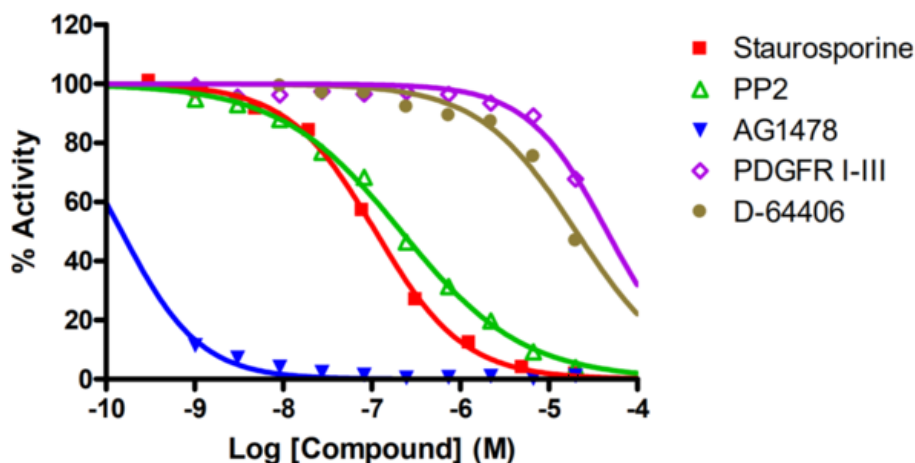
## Target Overview

|            |  |
|------------|--|
| CAT#       | EGFR (D770GY)  |
| Enzyme     | Human EGFR (D770GY)  |
| HGNC       | EGFR (D770GY)  |
| Synonym    | ERBB, mENA, ERBB1  |
| Substrates | Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml<br>ATP 10 μM |

## Assay Properties

| Reaction                 | Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> P-Substrate + ADP   |           |           |               |     |     |     |        |      |             |        |         |        |
|--------------------------|--|-----------|-----------|---------------|-----|-----|-----|--------|------|-------------|--------|---------|--------|
| Readout                  | Enzymatic activity   |           |           |               |     |     |     |        |      |             |        |         |        |
| Reference Compound IC50s | <table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>110</td> </tr> <tr> <td>PP2</td> <td>216</td> </tr> <tr> <td>AG1478</td> <td>&lt;1.0</td> </tr> <tr> <td>PDGFR I-III</td> <td>45,000</td> </tr> <tr> <td>D-64406</td> <td>20,000</td> </tr> </tbody> </table> | Compounds | IC50 (nM) | Staurosporine | 110 | PP2 | 216 | AG1478 | <1.0 | PDGFR I-III | 45,000 | D-64406 | 20,000 |
| Compounds                | IC50 (nM)  |           |           |               |     |     |     |        |      |             |        |         |        |
| Staurosporine            | 110  |           |           |               |     |     |     |        |      |             |        |         |        |
| PP2                      | 216  |           |           |               |     |     |     |        |      |             |        |         |        |
| AG1478                   | <1.0   |           |           |               |     |     |     |        |      |             |        |         |        |
| PDGFR I-III              | 45,000   |           |           |               |     |     |     |        |      |             |        |         |        |
| D-64406                  | 20,000   |           |           |               |     |     |     |        |      |             |        |         |        |
| Scientific Information   | <a href="http://www.uniprot.org/uniprot/P00533">P00533</a> ( <a href="http://www.uniprot.org/uniprot/P00533">http://www.uniprot.org/uniprot/P00533</a> )   |           |           |               |     |     |     |        |      |             |        |         |        |
| Screening Location       | Malvern, USA   |           |           |               |     |     |     |        |      |             |        |         |        |

## Reference compound IC50 for EGFR (D770GY)



Let's Discover Together.

© Aug 2022 **Reaction Biology**. All rights reserved.

[https://www.reactionbiology.com/datasheet/egfr\\_d770gy\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/egfr_d770gy_kin_malvern)

