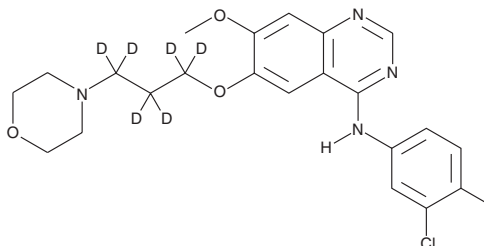


# PRODUCT INFORMATION



## Gefitinib-d<sub>6</sub> Item No. 22364

**CAS Registry No.:** 1228664-49-0  
**Formal Name:** N-(3-chloro-4-fluorophenyl)-7-methoxy-6-[3-(4-morpholinyl)propoxy-1,1,2,2,3,3-d<sub>6</sub>]-4-quinazolinamine  
**MF:** C<sub>22</sub>H<sub>18</sub>ClD<sub>6</sub>FN<sub>4</sub>O<sub>3</sub>  
**FW:** 452.9  
**Chemical Purity:** ≥95% (Gefitinib)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>6</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Gefitinib-d<sub>6</sub> is intended for use as an internal standard for the quantification of gefitinib (Item No. 13166) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Gefitinib-d<sub>6</sub> is supplied as a solid. A stock solution may be made by dissolving the gefitinib-d<sub>6</sub> in the solvent of choice, which should be purged with an inert gas. Gefitinib-d<sub>6</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of gefitinib-d<sub>6</sub> in ethanol is approximately 0.3 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Gefitinib-d<sub>6</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, gefitinib-d<sub>6</sub> should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Gefitinib-d<sub>6</sub> has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Gefitinib is an EGFR inhibitor (IC<sub>50</sub>s = 0.023-0.079 μM).<sup>1</sup> It inhibits colony formation of GEO colon, ZR-75-1 and MCF-10A breast, and OVCAR-3 ovarian cancer cell lines in soft agar assays (IC<sub>50</sub>s = 0.2-0.4 μM).<sup>2</sup> Gefitinib (0.1 to 1 μM) induces apoptosis and inhibits EGFR autophosphorylation in the same cells. *In vivo*, gefitinib (1.25, 2.5, and 5 mg/kg) reduces tumor volume and increases survival in a GEO mouse xenograft model. Formulations containing gefitinib have been used in the treatment of non-small cell lung cancer (NSCLC).

### References

1. Mendelsohn, J. and Baselga, J. The EGF receptor family as targets for cancer therapy. *Oncogene* **19**(56), 6550-6565 (2000).
2. Ciardiello, F., Caputo, R., Bianco, R., *et al.* Antitumor effect and potentiation of cytotoxic drugs activity in human cancer cells by ZD-1839 (Iressa™), an epidermal growth factor receptor-selective tyrosine kinase inhibitor. *Clin. Cancer Res.* **6**(5), 2053-2063 (2000).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/22/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM