

Product Information

Description

4-(4-(4-(3-fluorobenzoyloxy)-3-chlorophenylamino)quinazolin-6-yl)phenol, is a derivative of the quinazoline containing drug Lapatinib. Lapatinib reversibly blocks phosphorylation of the epidermal growth factor receptor (EGFR), ErbB2, and the Erk-1 and-2 and AKT kinases; it also inhibits cyclin D protein levels in human tumor cell lines and xenografts.

Molecular Formula:

C₂₇H₁₉ClFN₃O₂

Supplied As:

Solid

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

471.9 Da

Cas No.

1432450-51-5

Chemical Structure

