



ClariCELL™ FGFR1 Kinase Assay Service

4-Step Assay Validation

FGFR1 Expression in Cells

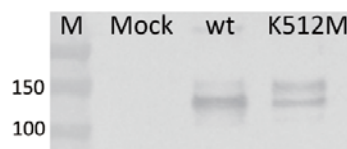


Figure 1: Wild type (wt) or kinase dead (K512M) FGFR1 was expressed transiently in 293 cells. Following cell lysis, an IP Western was performed with appropriate antibodies to capture and detect total FGFR1 protein.

FGFR1 Autophosphorylation in Cells

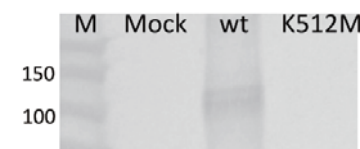


Figure 2: Wild type (wt) or kinase dead (K512M) FGFR1 was expressed transiently in 293 cells. An IP Western was performed with appropriate antibodies to capture and detect phospho-FGFR1 protein.

説明



ClariCELL™ FGFR1 Kinase Assay では、ヒト細胞内に transient に強制発現させた FGFR1 の自己リン酸化レベルを指標に、細胞内の FGFR1 に対する低分子化合物の阻害活性を検出します。

概要

HEK 293 細胞にヒト全長 FGFR1 を transient に導入します (アミノ酸配列は確認済みです)。FGFR1 の阻害剤をこの細胞に添加し、一定時間培養後、細胞を溶解させます。細胞内の FGFR1 をアッセイプレートの上に捕捉し、その自己リン酸化レベルを特異的の抗体を用いた ELISA によって定量します。ネガティブコントロールにはキナーゼ活性の無い FGFR1 [K512M] を導入した細胞中のリン酸化レベルの値を使用します。

Quantification of Phosphorylation

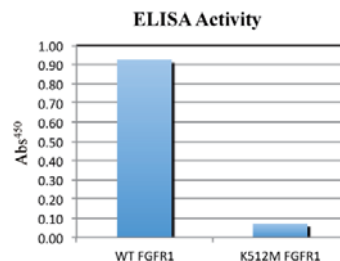


Figure 3: Wild type (wt) or kinase dead (K512M) FGFR1 was expressed transiently in 293 cells. Following cell lysis, an ELISA was performed to quantify the extent of autophosphorylation of FGFR1.

Reference Inhibitor Data

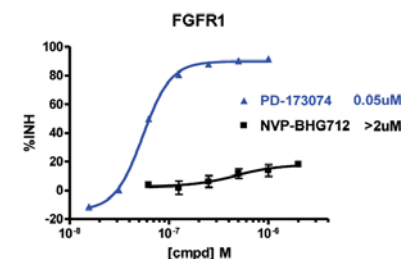


Figure 4: An autophosphorylation assay was performed in the presence of PD-173074, an FGFR1 inhibitor, and NVP-BHG712, a compound that is not expected to inhibit FGFR1. % inhibition data were plotted to determine EC₅₀ values.

サービスの詳細、最新情報はHPをご覧ください

カルナバイオサイエンス株式会社

〒650-0047 神戸市中央区港島南町1-5-5 BMA3F

TEL: 078-302-7091 (営業部直通) / FAX: 078-302-7086

E-mail: info@carnabio.com