

# PDGF receptor tyrosine kinase inhibitor IV

## Cell-permeable PDGFR inhibitor with antiproliferative properties.

A cell-permeable compound acting as an ATP-competitive and reversible inhibitor of PDGFR. PDGF receptor tyrosine kinase inhibitor IV displays potent antiproliferative properties in several human tumor cell lines.

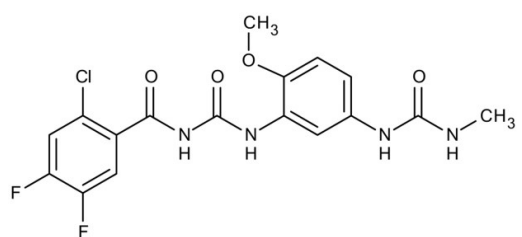
### Ordering Information

[Order Online »](#)

|                 |     |
|-----------------|-----|
| ENZ-CHM111-0005 | 5mg |
|-----------------|-----|

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                    |                                                                  |
|--------------------|------------------------------------------------------------------|
| Use/Stability      | As indicated on product label or CoA when stored as recommended. |
| Short Term Storage | -20°C                                                            |
| Long Term Storage  | -80°C                                                            |
| Shipping           | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                                          |
|------------------|------------------------------------------------------------------------------------------|
| Alternative Name | 3-Fluoro-N-(6,7-dimethoxy-2,4-dihydroindeno[1,2-c]pyrazol-3-yl)phenylamine, JNJ-10198409 |
| Appearance       | White powder solid.                                                                      |
| CAS              | 627518-40-5                                                                              |
| Couple Target    | PDGF receptor                                                                            |
| Couple Type      | Inhibitor                                                                                |
| Formula          | $C_{18}H_{16}FN_3O_2$                                                                    |
| MW               | 325.3                                                                                    |
| Purity           | ≥99% (HPLC)                                                                              |
| Solubility       | Soluble in DMSO.                                                                         |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)