

## Axitinib

**Catalog #:** 27064

**Lot #:** 220722

**Size:** 200 mg

**Structure:**

**CAS Registry #:** 319460-85-0

**Purity:** ≥99%

**Chemical Formula:** C<sub>22</sub>H<sub>18</sub>N<sub>4</sub>OS

**Molecular Weight:** 386.5 Da



**Description:** Axitinib (AG-013736) is an oral, potent, and selective inhibitor of vascular endothelial growth factor receptors 1 (VEGFR-1), 2 (VEGFR-2), and 3 (VEGFR-3), platelet derived growth factor receptor (PDGFR), and cKIT (CD117).

**Appearance:** White to off-white solid

**Solubility:** Soluble in DMSO at 33 mg/ml; soluble in ethanol at 1.7 mg/ml with warming; very poorly soluble in water; maximum solubility in plain water is estimated to be about 10-50 µM.

**Storage/Stability:** Store at or below –20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

### Reference:

1. Mancuso, M.R., et al. J. Clin. Invest. 116: 2610-2621 (2006).