

Vascular APIs

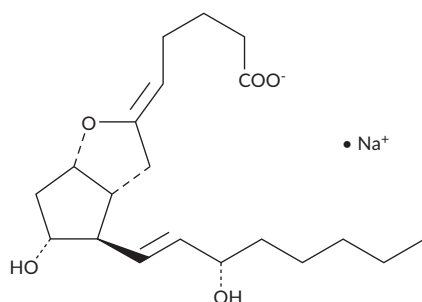
Each prostaglandin (PG) evokes distinct physiological effects through specific cell surface receptors. PGI₂ (prostacyclin) promotes smooth muscle relaxation and inhibits platelet activation through the IP receptor. Agonists of IP, including epoprostenol, iloprost, and treprostinil, are effective in relaxing pulmonary arterial smooth muscle and are used in the treatment of pulmonary hypertension.

- US FDA and EMA compliant
- Analytical standards of impurities and degradation products available



Commercial API Production

CGMP Epoprostenol sodium salt



CAS Number 61849-14-7

Availability GMP material is available, and the DMF is on file with the US FDA, Canada, Japan, Australia, and many EU member states.