

Animal Health APIs

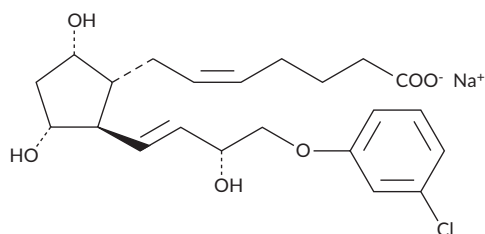
A stable and potent analog of prostaglandin $F_{2\alpha}$ ($PGF_{2\alpha}$), cloprostenol is used in the care of domestic mammals to induce luteolysis, initiate estrus, and synchronize estrus cycles. Both the racemic and active isomer are also used to induce parturition, terminate abnormal and normal pregnancies, and treat chronic endometriosis and ovarian luteal cysts.

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



Commercial API Production

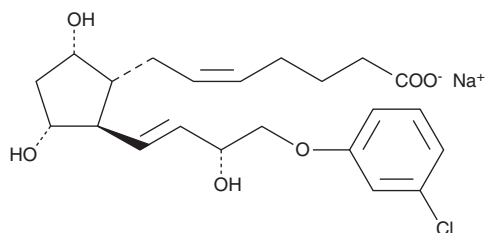
CGMP (±)-Cloprostenol sodium salt



CAS Number 55028-72-3

Availability GMP material, which conforms to the USP and BVP Monographs, is available. The DMF is on file with the US FDA, Canada, Russia, in almost all EU member states, and many other countries.

CGMP (+)-Cloprostenol sodium salt



CAS Number 62561-03-9

Availability GMP material is available, and the ASMF for (+)-Cloprostenol sodium salt is on file in the UK, Switzerland, and most EU countries.