

# Erythropoietin $\alpha$ (human), (recombinant)

## Recombinant human EPO that stimulates red blood cell production and increases oxygen capacity.

Erythropoietin (EPO) is a 34 kDa glycoprotein hormone produced by the kidney that promotes the formation of red blood cells in the bone marrow, thereby increasing the oxygen carrying capacity of blood in times of oxygen demand.

### Ordering Information

[Order Online »](#)

ADI-908-050-500	500U
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)

