

WesternSure® Pen Quick Reference Card

The WesternSure Pen is used to annotate visible protein ladders prior to chemiluminescent Western blot detection. The pen is optimized for LI-COR Imagers that detect chemiluminescence, including the C-DiGit® Blot Scanner, the Odyssey® Fc Imager, the Odyssey XF Imager, the Odyssey M Imager, and is suitable for use with film or other imaging systems. This unique marker delivers an ink which emits light when incubated with commonly used chemiluminescent substrates, including WesternSure PREMIUM. The ink is faintly visible for easy identification of marked membranes.

Ordering Information

WesternSure Pen [LI-COR P/N 926-91000]

www.licor.com/bio

Instructions

- Lightly touching the pen to the membrane should be enough to transfer ink to the membrane.
- Do not push down on the nib so hard that it creates an uneven surface on the membrane.
- Membranes may be annotated when damp after transfer, or when dry.
- Annotated membranes may be stored dry at ambient temperature or 4 °C for up to 1 week before starting the Western blot detection process.
- If ink is not flowing smoothly onto a damp membrane, trace over the band until it is annotated to the desired effect.

WesternSure™ Pen Quick Reference Card

Instructions

Please read all steps below prior to use.

1. Remove the pen cap just before use.

Caution: the pen will dry out if it is left uncapped or it is not sealed tightly before storage.

2. Test the pen on a piece of paper before writing on the blot to ensure that the pen is writing smoothly.
3. For best results, wick off excess buffer prior to annotation.

4. The pen can be used to annotate the visible ladder on the membrane at any point during the Western blot detection process **prior to the addition of substrate**; add other notations, as desired.
5. Upon addition of substrate containing luminol, the annotations will become visible when imaged or exposed to film.
6. If the blot is stripped and re-probed, annotate the visible ladder again, as outlined above.