

Urinary Indican Test Kit

Catalog Number: I1000N

DESCRIPTION

INDICAN (indoxyl sulfate) is a putrefaction product resulting from bacterial deconjugation of dietary tryptophan to indole in the small intestine. Indican levels are directly associated with bacterial activity in the intestines. Elevated levels indicate intestinal toxemia or overgrowth of anaerobic bacteria, putrefaction of undigested food in the bowels, stomach disorder (constipation, malabsorption), intestinal disorders and pancreatic insufficiency.

BioAssay Systems' Urinary Indican test kit is a less hazardous modification of the original Obermeyer indican reagent. It provides an accurate, simple, fast (15 min) and inexpensive screening tool for the identification of protein putrefaction in the gastrointestinal tract.

Precautions: reagents are for research use only. Not intended for diagnostic use. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

Product Information

Catalog No: I1000N

Number of Tests: 20 per package.

Contents:

- 20 Vials of Reagent
- 20 Droppers

Shipping/Storage: The kit is shipped and stored at room temperature. Protect from sunlight.

Expiry: see expiry date on the vial label.

TEST PROCEDURE

Caution: Corrosive! The Indican Reagent contains hydrochloric acid. When performing the test, wear gloves, goggles and protective clothing. In case of accidental contact with the reagent, immediately cover with baking soda and wash with ample amount of warm water.

Samples: No alcohol the night before; preferably a high protein meal the night before; no iodine or bile supplements taken in high doses 3 - 4 days prior to testing. Obtain first morning urine sample in a clean container. Urine Indican.

1. Unscrew the Cap off the Indican vial.
2. Using the dropper, carefully transfer 2 x 1mL of urine (each time fill dropper to the top 1-mL line) to the vial. Avoid spillage.
3. Replace cap, securely close the vial. Immediately invert the vial five times at 0, 5, 10 and 15 min.
4. Compare the color in the top solvent layer with the Indican Color Chart below. Record results and dispose the materials properly.

Expected Results

Clear or Light Yellow	0: Normal
Greenish Blue	1: Low Positive
Light Blue	2: Medium Positive
Blue	3: High Positive
Dark Blue	4: Very High Positive.

Indican Color Chart (Cat#: I1000N)

