

# Copper Gleam PCM or PC in acid copper baths

## Keywords

Brightener, acid copper plating, printed circuit boards, ion chromatography, IC, Dionex ICS-5000<sup>+</sup>

## Performance

The major active component in Copper Gleam™ PCM or Copper Gleam™ PC can be determined down to 0.01% in acid copper baths. The manufacturer's suggested concentration range is 0.4% to 1.0%.

## Application areas

- Acid copper plating
- Printed circuit boards

## Equipment

Thermo Scientific™ Dionex™ DX-500 Ion Chromatography system\*

\* Equivalent or improved results can be achieved using the Thermo Scientific™ Dionex™ ICS-5000<sup>+</sup> HPIC™ system with a Thermo Scientific™ Dionex™ ICS-Series VWD Variable Wavelength Detector.

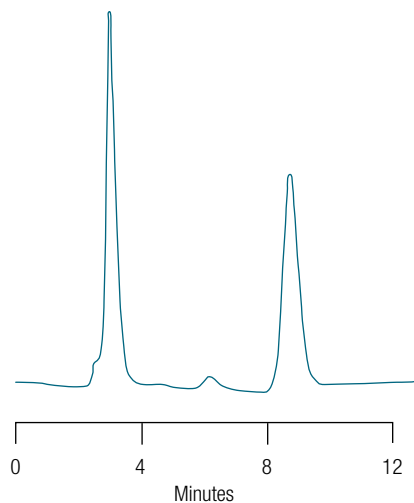


Figure 1. Pure Copper Gleam PCM at 254 nm.

### Conditions

Column:	Thermo Scientific™ Dionex™ IonPac™ NS1
Eluent:	10 mM H <sub>2</sub> SO <sub>4</sub> 10% CH <sub>3</sub> CN
Flow Rate:	1 mL/min
Sample Volume:	50 µL
Detection:	UV/Vis, 254 nm

### Sample preparation

Samples are injected directly using a 50 µL sample size.

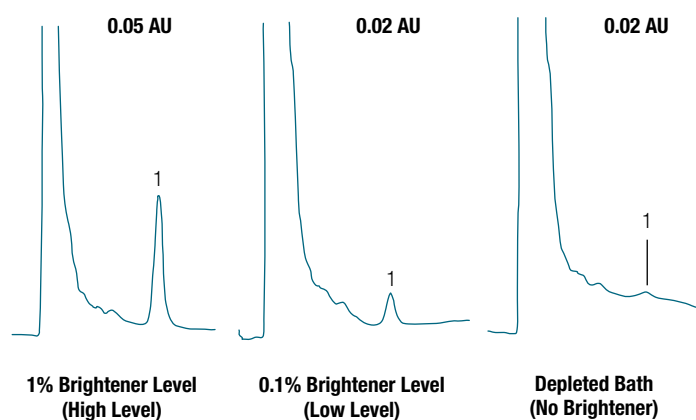


Figure 2. Determination of brightener in acid copper baths by ion chromatography.

Find out more at [thermofisher.com/IC](https://thermofisher.com/IC)

**ThermoFisher**  
SCIENTIFIC