

Product Information

CellLytic[™] P

Plant Cell Lysis/Extraction Reagent

Product Number **C 2360**

TECHNICAL BULLETIN

Product Description

The CellLytic[™] P reagent offers a convenient method for efficient plant cell lysis and protein solubilization, while avoiding protein degradation and interference with protein immunoreactivity and biological activity. CellLytic P reagent is a detergent-based reagent that may be used for the extraction of proteins from less than one gram of fresh or frozen leaves up to hundreds of grams, employing a short procedure (20 to 30 minutes). The CellLytic P reagent has been tested for use with tobacco, tomato, spinach and arabidopsis.

The resulting lysate is suitable for:

- Western blot analysis
- Protein staining such as Coomassie and silver staining
- Enzymatic assays: e.g. Kinase assays (tested for tyrosine and serine kinase activity) and phosphatase assays (tested for general and alkaline phosphatase)
- DNA-protein interaction assays (Gel-Shift)

Reagent

CellLytic P reagent is supplied ready-to-use. Each ml of reagent is sufficient for extraction of 0.5 to 1 gram of plant leaves.

Reagents and Equipment Required For Extraction but Not Provided

(Sigma product numbers are given where available)

- Protease Inhibitor Cocktail for Plants (P 9599)
- Test tubes
- Centrifuge or Microcentrifuge (Eppendorf 5417R or equivalent)
- Liquid nitrogen
- Mortar and pestle or homogenizer

Precautions and Disclaimer

Sigma's CellLytic P reagent is for laboratory use only. Not for drug, household or other uses.

Storage/Stability

Store at room temperature.

Procedure

To avoid protein degradation it is highly recommended to add Protease Inhibitor Cocktail to the CellLytic P reagent (diluted 1:100).

1. Grind fresh or frozen leaves to a fine powder with liquid nitrogen, using mortar and pestle: Add 2 ml of CellLytic P reagent to each gram of leaves and grind thoroughly with the pestle. Alternatively: Grind the leaves in blender or homogenizer together with the CellLytic P reagent. **Note:** For a higher concentration of extracted proteins, reduce the volume of CellLytic P reagent per gram of leaves.
2. Transfer the suspension to a tube(s).
3. Centrifuge for 10 minutes at 12,000 x g to pellet the cellular debris.
4. Transfer the supernatant to a new tube. At this step, the extract is suitable for immediate use. Alternatively, the extract may be stored at -70 °C. To improve protein stability, glycerol may be added a final concentration of 10 to 25 % prior to freezing as long as this will not interfere with your application.

YA/EB/AC 8/02

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.