

DNAbsolute

Column-free DNA extraction kit



DNAbsolute is an ionic liquid that immediately binds and precipitates DNA strands when added to your lysate. It provides a simple method combining safety, speed and sensitivity.

Key features

1. Safe & simple

No need for any columns or harmful chemicals, highly amenable to automated workflows & on-field applications

2. Versatile

Extract DNA from a variety of samples, including some of the most challenging ones (i.e. insects, plants & corals)

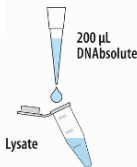


3. Extreme sensitivity

Efficient DNA recovery from reduced starting materials, including individual specimens and even down to a single leg!

How does it work?

1. Lyse

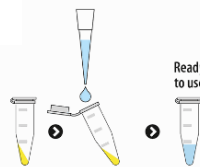


Centrifuge
5 min

3. Wash & resuspend



Centrifuge
5 min



2. Add DNAbsolute

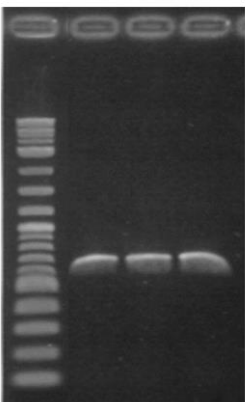
Kit contents:

- DNAbsolute reagent, 3M
- PBS 150 mM, pH 7.2
- Tris 25 mM, pH 8
- 2x Lysis buffers

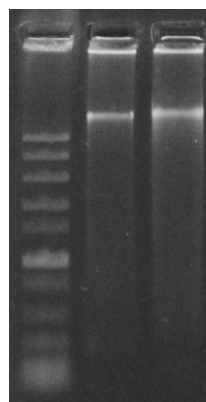


Results

Kit DNAbs. DNAbs.



Extraction of PCR-grade DNA from *Drosophila* using DNAbsolute



Extraction of highly pure DNA from coral specimens using DNAbsolute

Credits: Gabriel Johnson - Smithsonian Institution, Museum Support Center, Maryland, US