

Technical Data Sheet

**iD PAGE GELS**  
**ID-PAYYYY-XXX**

Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics.

**Products**

Cat #	Description	Size
ID-PA4201-010	iD PAGE Gel, 4-20%, 10 wells	1 Kit – 10 gels
ID-PA4201-012	iD PAGE Gel, 4-20%, 12 wells	
ID-PA4201-015	iD PAGE Gel, 4-20%, 15 wells	
ID-PA4121-010	iD PAGE Gel, 4-12%, 10 wells	
ID-PA4121-012	iD PAGE Gel, 4-12%, 12 wells	
ID-PA4121-015	iD PAGE Gel, 4-12%, 15 wells	
ID-PA8161-010	iD PAGE Gel, 8-16%, 10 wells	
ID-PA8161-012	iD PAGE Gel, 8-16%, 12 wells	
ID-PA8161-015	iD PAGE Gel, 8-16%, 15 wells	
ID-PA0121-010	iD PAGE Gel, 12%, 10 wells	
ID-PA0121-012	iD PAGE Gel, 12%, 12 wells	
ID-PA0121-015	iD PAGE Gel, 12%, 15 wells	
ID-PA0101-010	iD PAGE Gel, 10%, 10 wells	
ID-PA0101-012	iD PAGE Gel, 10%, 12 wells	
ID-PA0101-015	iD PAGE Gel, 10%, 15 wells	
ID-PA0081-010	iD PAGE Gel, 8%, 10 wells	
ID-PA0081-012	iD PAGE Gel, 8%, 12 wells	
ID-PA0081-015	iD PAGE Gel, 8%, 15 wells	

**Description**

The iD PAGE Gels are 1mm thick precast mini polyacrylamide gels perfectly suited for large loading volumes (up to 80 µl). They are casted in a neutral pH buffer that minimizes the hydrolysis of polyacrylamide and results in extra gel stability.

Because they are manufactured without SDS, the iD PAGE Gels are ideal for both SDS-PAGE and native electrophoresis depending on the running buffer and transfer buffer used.

The design of the cassette was optimised to provide a better band resolution and a significant improvement of the sample distribution in the loading wells.

The proprietary gel-casting techniques provide an excellent batch-to-batch consistency and guarantee a reliable migration pattern. Using specially formulated Tris-MOPS running buffer, these gels enable proteins to be separated quickly and easily for subsequent detection by staining or western blotting.

**Compatibility**

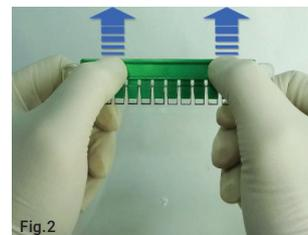
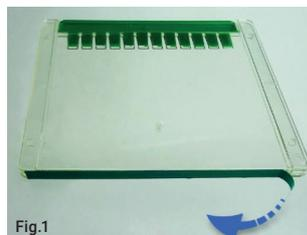
The iD PAGE Gels are compatible with the following Gel Tanks:

- Bio-Rad Mini-PROTEAN® II & 3
- Bio-Rad Mini-PROTEAN® Tetra System
- LONZA PAGE® Minigel Chamber
- Hoefer Mighty Small (SE 260/SE 250)
- Hoefer Tall Mighty Small (SE 280)
- Invitrogen Novex XCell I, II, & Surelock® (Use with Eurogentec Gel Tank Adapter Plates)

**Storage**

Store at 4-25°C. Gels are stable for up to 18 months if stored at 2-8°C

**Instruction for use**



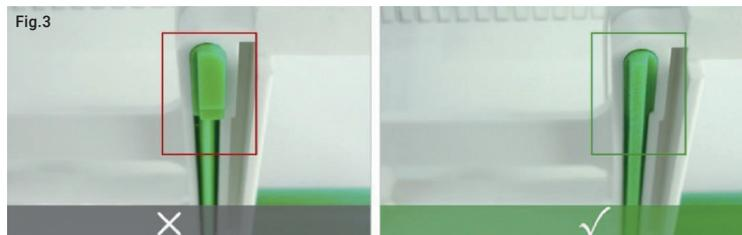
1) Remove the iD PAGE Gel from the package, peel the sealing tape at the bottom of the gel cassette. (Fig 1.)

2) Remove the comb from the gel cassette gently (Fig 2.)

3) Insert the gel into the gel running apparatus. Please refer to the apparatus manufacturer's instructions.

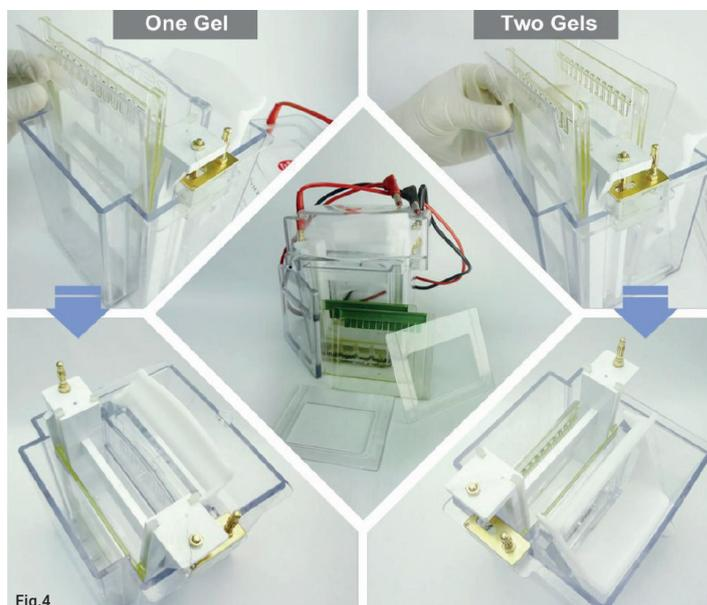
**NOTE :**

**Using Bio-Rad Mini-PROTEAN® Tetra System:** remove the gasket from the inner frame, turn it around so the flat side is facing outwards and insert the gasket back into the inner frame (Fig 3).



**Using Invitrogen Novex Mini-Cell tanks:** Adapters provided in the package are needed since the iD PAGE Gel cassette is thinner than the Invitrogen NuPAGE® gel cassette.

**Using Invitrogen Novex® Mini-Cell:** Please refer to figure 4 (Fig 4).



## Technical Data Sheet

4) Pour sufficient 1x iD MOPS running buffer into the inner tank of the gel running apparatus to cover the sample wells by 5-7 mm. Fill the outer tank with the same running buffer to ensure proper cooling. For best results, the buffer in the outer tank should be above the top level of the sample wells.

*NOTE: Do NOT use tris-glycine running buffer with the iD PAGE Gels.*

5) Rinse the sample wells thoroughly with 1x running buffer to remove air bubbles and displace any storage buffer before loading the samples.

6) Prepare your samples

*FOR NATIVE PAGE, please note that the iD PAGE Gels are precast without SDS which is conducive for native PAGE. Protein samples should be prepared in a non-reducing, non-denaturing sample buffer, to maintain the proteins' secondary structure and native charge. The mobility of the protein depends on the size and shape of the protein as well as its net charge.*

7) Load your samples

Make sure the loading tip is vertically inserted into the loading well for optimal results. Protein overloading will cause smearing and distortion. Excessive loading of proteins with free carbohydrates may also lead to band distortion or failure of the protein to penetrate the gel.

8) Run the gel (140 V; 40-60 minutes depending on the protein size).

9) After protein migration, remove the gel from the gel tank and open the cassette carefully by inserting the cassette opener into the gap between the two plates. A cracking sound may be heard as you open the cassette. Please wear protective goggles to avoid eye contacts or damages (Fig 5).

10) Loosen the gel from the plate with water and gently remove. Please dispose used cassettes as non-hazardous medical waste.

### Staining

All standard SDS staining procedures can be used with the iD PAGE gels. When using commercially available staining reagents and device, follow the manufacturer's instructions.

### Protein Transfer

All standard transferring procedures can be used with the iD PAGE Gels.

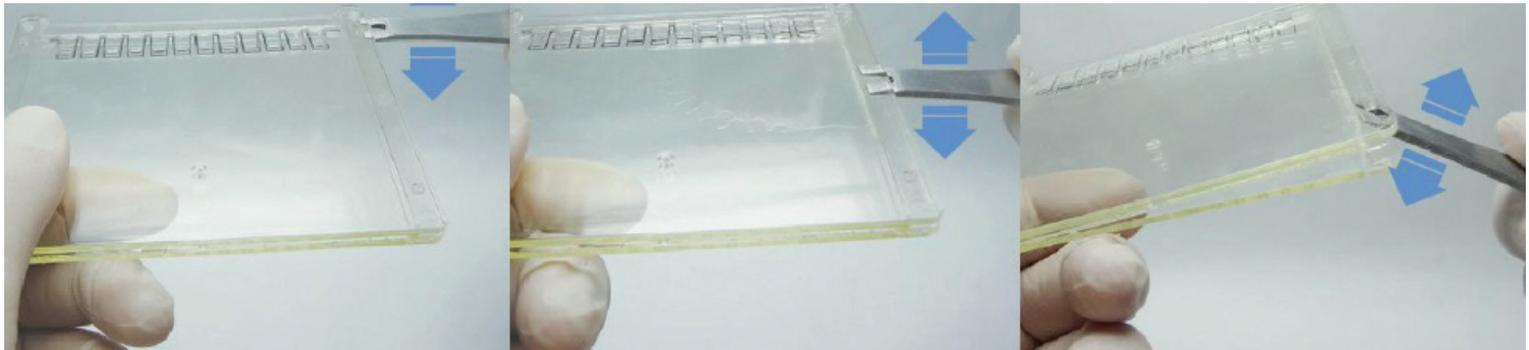


Fig.5

**For further information please contact our Customer Help Desk:**

**For Europe:**

E-mail: info@eurogentec.com  
Tel: +32 4 372 76 65

**For USA:**

E-mail: service@anaspec.com  
Tel: +1 510 791 9560