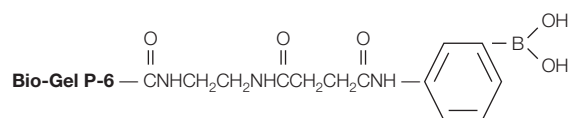


Affi-Gel® Boronate Affinity Gel to Separate Ribonucleotides, Ribonucleosides, Sugars, Catecholamines, and Coenzymes

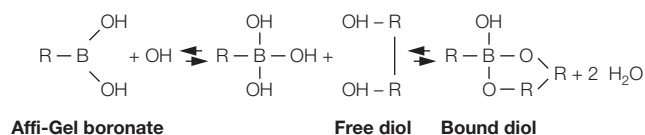
Introduction

Low molecular weight molecules such as nucleotides, nucleosides, catecholamines, coenzymes, and sugars can be separated very efficiently by chromatography on Affi-Gel boronate gel. Affi-Gel boronate gel is a boronate-derivatized polyacrylamide gel with an affinity for adjacent *cis* hydroxyl (*cis*-diol) groups (Figure 1). It effectively separates molecules, such as AMP and cyclic AMP (cAMP), which differ only in the presence or absence of such a group. It also separates *cis*-diol-containing molecules such as cytosine, uridine, and adenosine from one another — all are bound, but their differing affinities permit separate elution (Figure 2A). Bio-Gel® P-6 beads, the gel beads derivatized to make Affi-Gel boronate gel, have a molecular weight exclusion limit of 6,000. The binding capacity of Affi-Gel boronate gel is, therefore, extremely high for low molecular weight compounds.

The development and use of boronate affinity media stems from the observation that boronate forms complexes with sugars possessing coplanar *cis*-diol groups (Boeseken 1949). The structure of the boronate affinity group in Affi-Gel boronate gel is:

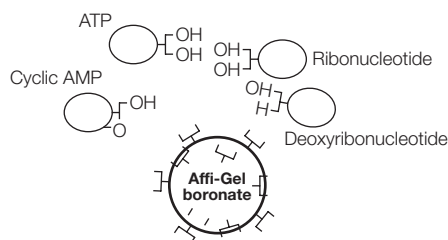


The mechanism of binding has been suggested to be reversible formation of a five-member ring (Lorand et al. 1959).



Boronate derivatives on various solid supports have been investigated and applied (Table 1). These derivatives have been variously termed dihydroxy boryl, boric acid gel, and boronic acid gel, all referring to *m*-aminophenyl boronic acid immobilized to solid supports. Supports have included aminoethyl cellulose, porous glass, methacrylate gel, and Bio-Gel P polyacrylamide gel.

Sample application



Retention of *cis*-diol compounds

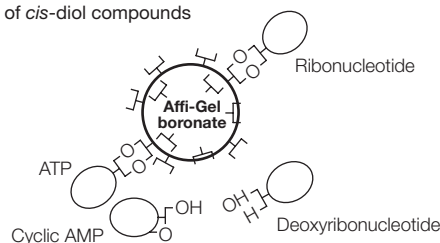


Fig. 1. Separation by affinity for *cis*-diols.

Affi-Gel boronate gel offers a variety of advantages as a support for the applications listed in Table 1:

- High boronate capacity — 1.05 ± 0.15 meq/g
- High capacity for diols — binds 130 μmol sorbitol/ml gel*
- Spherical polyacrylamide matrix — excellent flow rate, resolution, and reproducibility
- Low nonspecific absorption, optimum porosity — low exclusion limit precludes significant binding of high molecular weight molecules, while allowing rapid equilibration and binding of low and medium molecular weight molecules

Preparation and Use

The gel is supplied as dried material and must be hydrated in an appropriate buffer. Buffer choice is dependent upon application (Table 1). Buffers such as Tris should be avoided, since they can adversely affect binding capacity (Schott et al. 1973). The presence of Mg^{2+} may enhance binding. In general, binding is effected at $\text{pH} > 7.5$ and elution at $\text{pH} < 6.5$. Boric acid, sorbitol, or mannitol can also be used for elution.

* Dr R Hurlbert, personal communication.

Table 1. Applications using boronate affinity media.

Application	Application Buffer	Molecules in V ₀	Retained Molecules	Elution Buffer	Support Used	References
Adenylate cyclase assay	HEPES pH 7.5 + MgCl ₂	cAMP	ATP, AMP, and adenosine	0.05 M NaOAc	Bio-Gel P-150 boronate gel	Hageman and Kuehn 1977
Isolation of catecholamines from urine	0.1 M phosphate pH 7.0 + EDTA	Other urine components and DOPA	Norepinephrine, epinephrine	0.025 N HCl	Boric acid gel (Sigma-Aldrich)	Higa et al. 1977
Modified nucleosides in urine	0.25 M NH ₄ OAc, pH 8.8	Thymidine, adenine	Pseudouridine	0.1 M HCOOH	Bio-Gel P-2 gel, 200–400 mesh, boronate	Davis et al. 1977 Uziel et al. 1976
Separation of mono- and oligonucleotides	Triethyl ammonium	Deoxyribo-nucleotides	Ribonucleotides	H ₂ O and others	Dihydroxyboryl methacrylate	Schott et al. 1973
Sugars	0.05 M N-methyl-morpholinium-Cl, pH 7.5 + 1 M NaOAc	Erythritol, adonitol, sucrose, and D-glucose	L-arabinitol, xylitol, D-mannitol, dulcitol, sorbitol, and D-fructose*	Isocratic run, elution buffer same as application buffer	Dihydroxyboryl cellulose	Weith et al., 1970
Assay of ribonucleotide reductase	Tris, MgCl ₂	dUMP, dCMP	CDP	Citrate	Dihydroxyboryl cellulose Dowex 1(AG) resin	Moore et al. 1974

* In order of increasing retention by support.

Regeneration and Storage

Regeneration of the gel is facilitated by washing with starting buffer if a low-pH elution buffer has been used. If boric acid buffer or diol solutions have been used for elution, the gel must first be washed with 0.1 M acetic acid before reequilibrating with starting buffer.

Hydrated gel can be stored at 4°C under neutral pH conditions for up to 2 months; dry material has excellent long-term stability.

Typical Procedure

Stepwise elution is typical for separation of a ribonucleotide (such as GMP) from a cyclic nucleotide (cGMP) (Figure 2B):

1. Prepare a 1 x 4 cm column of Affi-Gel boronate in 0.1 M HEPES, pH 8.5.
2. Load a 200 µl solution containing 1 µmol each cGMP and GMP in the starting buffer onto the column.
3. Wash the column with 5 column volumes of starting buffer at a flow rate of 1 ml/min and collect the cyclic nucleotide.
4. Elute the ribonucleotide with 0.1 M NaPO₄, pH 6.0.

References

- Boeseken J, The use of boric acid for the detection of the configuration of carbohydrates, *Advan Carbohydr Chem* 4, 189–210 (1949)
- Davis GE et al., High-performance liquid chromatographic separation and quantitation of nucleosides in urine and some other biological fluids, *Clin Chem* 23, 1427–1435 (1977)
- Hageman JH and Kuehn GD, Assay of adenylate cyclase by use of polyacrylamide-boronate gel columns, *Anal Biochem* 80, 547–554 (1977)
- Higa S et al., Isolation of catecholamines in biological fluids by boric acid gel, *Anal Biochem* 77, 18–24 (1977)
- Lorand JP and Edwards JO, Polyol complexes and structure of the benzenboronate ion, *J Org Chem* 24, 769–774 (1959)
- Moore EC et al., Separation of ribonucleotides and deoxyribonucleotides on columns of borate covalently linked to cellulose. Application to the assay of ribonucleoside diphosphate reductase, *Biochemistry* 13, 2904–2907 (1974)

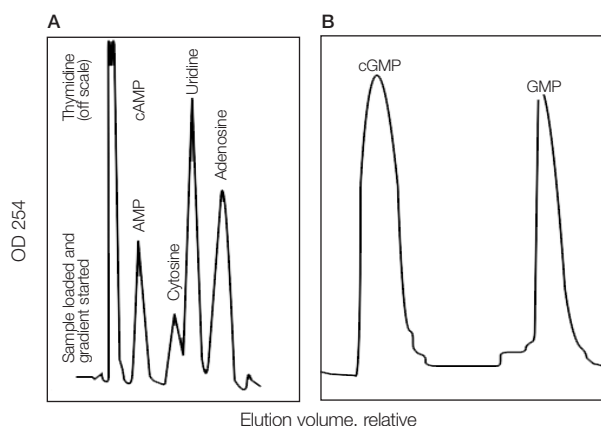


Fig. 2. Examples of separations using Affi-Gel boronate. A, separation of thymidine, cAMP, AMP, adenosine, cytosine, and uridine by a typical gradient elution. A sample containing 0.5 µmol of each molecule was loaded on an equilibrated 0.7 x 45 cm bed of Affi-Gel boronate gel; 0.1 M phosphate was applied after the first peak emerged. The flow rate was 40 ml/hr. B, separation of cGMP and GMP. A sample containing 1.0 µmol cGMP and GMP was loaded onto a 1 x 4 cm column equilibrated with 0.1 M HEPES, pH 8.5. GMP was eluted with 0.1 M NaPO₄, pH 6.0, and detected by a UV monitor.

Schott H et al., A dihydroxyboryl-substituted methacrylic polymer for the column chromatographic separation of mononucleotides, oligonucleotides, and transfer ribonucleic acid, *Biochemistry* 12, 932–938 (1973)

Uziel M et al., Modified nucleosides in urine: selective removal and analysis, *Clin Chem* 22, 1451–1455 (1976)

Weith HL et al., Synthesis of cellulose derivatives containing the dihydroxyboryl group and a study of their capacity to form specific complexes with sugars and nucleic acid components, *Biochemistry* 9, 4396 (1970)

We wish to thank Dr Robert Hurlbert, Biochemistry Department, MD Anderson Hospital, Houston, TX for his substantial contribution to the development of the chromatographic material described here.

Dowex is a trademark of the Dow Chemical Company.

Ordering Information

Catalog #	Description
153-6103	Affi-Gel Boronate Gel, 5 g
153-6104	Affi-Gel Boronate Gel, 50 g

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 2527 3454 **Canada** (905) 712-2771 **China** (86 21) 6426 0808 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84-0 **Greece** 30 210 777 4396 **Hong Kong** (852) 2789 3300 **Hungary** 36 1 455 8800 **India** (91-124)-2398112/3/4, 5018111, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 55-52-00-05-20 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 331 99 99 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-64153188 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 52 00 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** (886 2) 2578 7189/2578 7241 **United Kingdom** 020 8328 2000