



# Use of Aminex HPLC Columns for Carbohydrate Analysis

## Publications List

### 2007

**Chheda JN and Dumesic JA (2007).**

An overview of dehydration, aldol-condensation and hydrogenation processes for production of liquid alkanes from biomass-derived carbohydrates.

Catal Today 123, 59–70.

**Román-Leshkov Y et al. (2007).**

Production of dimethylfuran for liquid fuels from biomass-derived carbohydrates.

Nature 447, 982–985.

### 2010

**Arrizon J et al. (2010).**

Comparison of the water-soluble carbohydrate composition and fructan structures of *Agave tequilana* plants of different ages.

Food Chem 122, 123–130.

### 2011

**Weingarten R et al. (2011).**

Design of solid acid catalysts for aqueous-phase dehydration of carbohydrates:

The role of Lewis and Brønsted acid sites.

J Catal 279, 174–182.

## 2012

**Sluiter A et al. (2012).**

Determination of Structural Carbohydrates and Lignin in Biomass.

U.S. Department of Energy, Office of Energy Efficiency & Renewable Energy, National Renewable Energy Laboratory. Technical Report NREL/TP-510-42618.

**Yang Y et al. (2012).**

Conversion of carbohydrates and lignocellulosic biomass into 5-hydroxymethylfurfural using  $\text{AlCl}_3 \cdot 6\text{H}_2\text{O}$  catalyst in a biphasic solvent system.

Green Chem 14, 509–513.

## 2017

**An R et al. (2017).**

Efficient one-pot synthesis of n-butyl levulinate from carbohydrates catalyzed by  $\text{Fe}_2(\text{SO}_4)_3$ .

J Energy Chem 26, 556–563.

**Hafid HS et al. (2017).**

Over production of fermentable sugar for bioethanol production from carbohydrate-rich Malaysian food waster via sequential acid-enzymatic hydrolysis pretreatment.

Waste Manag 67, 95–105.

**Le C and Stuckey DC (2017).**

Impact of feed carbohydrates and nitrogen source on the production of soluble microbial products (SMPs) in anaerobic digestion.

Water Res 122, 10–16.

**Lin H et al. (2017).**

Conversion of carbohydrates into 5-hydroxymethylfurfural in a green reaction system of  $\text{CO}_2$ -water-isopropanol.

AIChE Journal 63, 257–265.

**Liu ZH et al. (2017).**

Synergistic maximization of the carbohydrate output and lignin processability by combinatorial pretreatment.

Green Chem 19, 4,939–4,955.

**Peng K et al. (2017).**

Hydrothermally stable Nb-SBA-15 catalysts applied in carbohydrate conversion to 5-hydroxymethyl furfural.

Mol Catal 441, 72–80.

**Wang S et al. (2017).**

Conversion of C5 carbohydrates into furfural catalyzed by a Lewis acidic ionic liquid in renewable  $\gamma$ -valerolactone.

Green Chem 19, 3,869–3,879.

## 2018

**Jiang L et al. (2018).**

Direct catalytic conversion of carbohydrates to methyl levulinate: Synergy of solid Brønsted acid and Lewis acid.  
Appl Catal B 220, 589–596.

**Li K et al. (2018).**

Quasi-homogeneous carbocatalysis for one-pot selective conversion of carbohydrates to 5-hydroxymethylfurfural using sulfonated graphene quantum dots.  
Carbon 136, 224–233.

**Vénica CI et al. (2018).**

Effect of the carbohydrates composition on physiochemical parameters and metabolic activity of starter culture in yogurts.  
LWT 94, 163–171.

**Yin W et al. (2018).**

Hydrotreatment of the carbohydrate-rich fraction of pyrolysis liquids using bimetallic Ni based catalyst: Catalyst activity and product property relations.  
Fuel Process Technol 169, 258–268.

## 2020

**Li X et al. (2020).**

Functional validation of two fungal subfamilies in carbohydrate esterase family 1 by biochemical characterization of esterases from uncharacterized branches.  
Front Bioeng Biotechnol 8, 694.

**Magri A et al. (2020).**

Production of fructooligosaccharides by *Bacillus subtilis natto* CCT7712 and their antiproliferative potential.  
J Appl Microbiol 128, 1,414–1,426.

**Viswanathan MB et al. (2020).**

Variability in structural carbohydrates, lipid composition, and cellulosic sugar production from industrial hemp varieties.  
Ind Crops Prod 157, 112,906.

Visit [bio-rad.com/AminexColumns](https://www.bio-rad.com/AminexColumns) for a wide range of videos, technical reports, published findings, and other resources describing Aminex Columns.

BIO-RAD and AMINEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.  
All trademarks used herein are the property of their respective owner.



**Bio-Rad  
Laboratories, Inc.**

---

Life Science  
Group

**Website** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

