



# Use of Aminex HPLC Columns in Biofuels Research

## Publications List

### 2010

**Pourbafrani M et al. (2010).**

Production of biofuels, limonene and pectin from citrus wastes.  
Bioresour Technol 101, 4,246–4,250.

**Qureshi N et al. (2010).**

Production of butanol (a biofuel) from agricultural residues: Part I — Use of barley straw hydrolysate.  
Biomass and Bioenergy 34, 559–565.

### 2013

**Kwon EE et al. (2013).**

Sequential co-production of biodiesel and bioethanol with spent coffee grounds.  
Bioresour Technol 136, 475–480.

### 2014

**Cai CM et al. (2014).**

Coupling metal halides with a co-solvent to produce furfural and 5-HMF at high yields directly from lignocellulosic biomass as an integrated biofuels strategy.  
Green Chem 16, 3,819–3,829.

### 2016

**Higgins BT et al. (2016).**

Cofactor symbiosis for enhanced algal growth, biofuel production, and wastewater treatment.  
Algal Research 17, 308–315.

## 2017

**Bhatia SK et al. (2017).**

Microbial biodiesel production from oil palm biomass hydrolysate using marine *Rhododoccus* sp. YHY01. *Bioresour Technol* 233, 99–109.

**Das L et al. (2017).**

Industrial hemp as a potential bioenergy crop in comparison with kenaf, switchgrass and biomass sorghum. *Bioresour Technol* 244, 641–649.

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

Direct conversion of biomass components to the biofuel methyl levulinate catalyzed by acid-base bifunctional zirconia-zeolites. *Applied Catalysis B: Environmental* 200, 182–191.

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

Hydrophobic Pd nanocatalysts for one-pot and high-yield production of liquid furanic biofuels at low temperatures. *Applied Catalysis B: Environmental* 215, 18–27.

**Maneechakr P and Karnjanakom S (2017).**

Catalytic transformation of furfural into bio-based succinic acid via ultrasonic oxidation using  $\beta$ -cyclodextrin-SO<sub>3</sub>H carbon catalyst: A liquid biofuel candidate. *Energy Conversion and Management* 154, 299–310.

**Ranganathan A et al. (2017).**

Utilizing anaerobic fungi for two-stage sugar extraction and biofuel production from lignocellulosic biomass. *Front Microbiol* 8, 635.

**Sun J et al. (2017).**

One-pot integrated biofuel production using low-cost biocompatible protic ionic liquids. *Green Chem* 19, 3,152–3,163.

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

Utilization of biodiesel by-product as substrate for high-production of  $\beta$ -farnesene via relatively balanced mevalonate pathway in *Escherichia coli*. *Bioresour Technol* 243, 228–236.

## 2018

**Calicioglu O and Brennan RA (2018).**

Sequential ethanol fermentation and anaerobic digestion increases bioenergy yields from duckweed. *Bioresour Technol* 257, 344–348.

**Farmanbordar S et al. (2018).**

Municipal solid waste as a suitable substrate for butanol production as an advanced biofuel. *Energy Conversion and Management* 157, 396–408.

**Mahmoodi P et al. (2018).**

Efficient conversion of municipal solid waste to biofuel by simultaneous dilute-acid hydrolysis of starch and pretreatment of lignocelluloses. *Energy Conversion and Management* 166, 569–578.

**Min D et al. (2018).**

Combination of hydrothermal pretreatment and sodium hydroxide post-treatment applied on wheat straw for enhancing its enzymatic hydrolysis. *Cellulose* 25, 1,197–1,206.

**Papadaki A et al. (2018).**

Fumaric acid production using renewable resources from biodiesel and cane sugar production processes. *Environ Sci Pollut Res Int* 25, 35,960–35,970.

**Sundstrom E et al. (2018).**

Demonstrating a separation-free process coupling ionic liquid pretreatment, saccharification, and fermentation with *Rhodospiridium toruloides* to produce advanced biofuels. *Green Chem* 20, 2,870–2,879.

## 2019

**Bhatia SK et al. (2019).**

Bioconversion of barley straw lignin into biodiesel using *Rhodococcus* sp. YHY01. *Bioresour Technol* 289, 121,704.

**Wang P et al. (2019).**

Towards comprehensive lignocellulosic biomass utilization for bioenergy production: Efficient biobutanol production from acetic acid pretreated switchgrass with *Clostridium saccharoperbutylacetonicum* N1-4. *Applied Energy* 236, 551–559.

**Wang Z et al. (2019).**

Improving ethanol yields with deacetylated and two-stage pretreated corn stover and sugarcane bagasse by blending commercial xylose-fermenting and wild type *Saccharomyces* yeast. *Bioresour Technol* 282, 103–109.

## 2020

### Dong L et al. (2020).

High-level expression of highly active and thermostable trehalase from *Myceliophthora thermophila* in *Aspergillus niger* by using the CRISPR/Cas9 tool and its application in ethanol fermentation. *J Ind Microbiol Biotechnol* 47, 133–144.

### Ebrahimian F and Karimi K (2020).

Efficient biohydrogen and advanced biofuel coproduction from municipal solid waste through a clean process. *Bioresour Technol* 300, 122,656.

### Hosseini A et al. (2020).

Efficient superantioxidant and biofuel production from microalga *Haematococcus pluvialis* via a biorefinery approach. *Bioresour Technol* 306, 123,100.

### Patel AK et al. (2020).

A sustainable mixotrophic microalgae cultivation from dairy wastes for carbon credit, bioremediation and lucrative biofuels. *Bioresour Technol* 313, 123,681.

### Wang B et al. (2020).

Bioenergy recovery from wastewater accelerated by solar power: Intermittent electro-driving regulation and capacitive storage biomass. *Water Res* 175, 115,696.

### Wang X et al. (2020).

Sustainable and stepwise waste-based utilisation strategy for the production of biomass and biofuels by engineered microalgae. *Environ Pollut* 265, 114,854.

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](https://www.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

